

THE COMPANIES ACTS 1985 to 2006

**ARTICLES OF ASSOCIATION
OF
CYPROTEX PLC**

Company Number 4311107

Amended by Special Resolution passed on 14 July 2009

CONTENTS

| Article No | Page |
|--|-------------|
| PRELIMINARY | 1 |
| 1 No other regulations to apply..... | 1 |
| 2 Interpretation | 1 |
| 3 Form of resolution | 4 |
| 4 Registered office | 4 |
| SHARE CAPITAL | 5 |
| 5 Authorised share capital..... | 5 |
| 6 Uncertificated shares..... | 5 |
| 7 Allotment..... | 8 |
| 8 Redeemable shares | 8 |
| 9 Power to attach rights..... | 8 |
| 10 Share warrants to bearer..... | 8 |
| 11 Commission and brokerage..... | 9 |
| 12 Trusts not to be recognised | 9 |
| ALTERATION OF SHARE CAPITAL | 9 |
| 13 Increase, consolidation, cancellation and sub-division..... | 9 |
| 14 Fractions of shares..... | 10 |
| 15 Reduction of capital..... | 11 |
| 16 Purchase of own shares | 11 |
| VARIATION OF CLASS RIGHTS | 11 |
| 17 Sanction to variation..... | 11 |
| 18 Class meetings..... | 11 |
| 19 Deemed variation | 12 |
| EVIDENCE OF TITLE | 12 |
| 20 Right to share certificates..... | 12 |
| 21 Replacement certificates | 13 |

| | |
|---|----|
| LIEN ON SHARES | 13 |
| 22 Lien on shares not fully paid..... | 13 |
| 23 Enforcement of lien by sale | 13 |
| 24 Application of proceeds of sale | 14 |
| CALLS ON SHARES | 14 |
| 25 Calls | 14 |
| 26 Liability of joint holders | 15 |
| 27 Interest on calls | 15 |
| 28 Rights of member when call unpaid | 15 |
| 29 Sums due on allotment treated as calls | 15 |
| 30 Power to differentiate..... | 15 |
| 31 Payment in advance of calls | 15 |
| 32 Delegation of power to make calls | 16 |
| FORFEITURE OF SHARES | 16 |
| 33 Notice if call not paid | 16 |
| 34 Forfeiture for non-compliance | 16 |
| 35 Notice after forfeiture | 16 |
| 36 Forfeiture may be annulled | 17 |
| 37 Surrender | 17 |
| 38 Disposal of forfeited shares | 17 |
| 39 Effect of forfeiture | 17 |
| 40 Extinction of claims..... | 18 |
| 41 Evidence of forfeiture..... | 18 |
| TRANSFER OF SHARES | 18 |
| 42 Transfer of certificated shares..... | 18 |
| 43 Transfer of uncertificated shares..... | 18 |
| 44 Right to refuse registration..... | 19 |
| 45 Notice of refusal..... | 19 |

| | | |
|----|--|----|
| 46 | Closing of register | 20 |
| 47 | Fees on registration..... | 20 |
| 48 | Transfers by renunciation | 20 |
| | TRANSMISSION OF SHARES..... | 20 |
| 49 | On death..... | 20 |
| 50 | Election of person entitled by transmission | 20 |
| 51 | Rights on transmission | 21 |
| | DESTRUCTION OF DOCUMENTS..... | 21 |
| 52 | Destruction of documents..... | 21 |
| | GENERAL MEETINGS..... | 22 |
| 53 | Annual general meetings..... | 22 |
| 54 | Convening of general meetings | 22 |
| 55 | Notice of general meetings | 22 |
| 56 | Omission to send notice | 24 |
| | PROCEEDINGS AT GENERAL MEETINGS | 24 |
| 57 | Quorum | 24 |
| 58 | If quorum not present | 24 |
| 59 | Chairman | 25 |
| 60 | Directors may attend and speak | 25 |
| 61 | Power to adjourn | 25 |
| 62 | Notice of adjourned meeting..... | 25 |
| 63 | Business of adjourned meeting..... | 26 |
| 64 | Accommodation of members and security arrangements..... | 26 |
| | VOTING | 27 |
| 65 | Method of voting..... | 27 |
| 66 | Chairman's declaration conclusive on show of hands | 27 |
| 67 | Objection to error in voting..... | 27 |
| 68 | Amendment to resolutions..... | 28 |

| | | |
|----|--|----|
| 69 | Procedure on a poll | 28 |
| 70 | Votes of members | 29 |
| 71 | Failure to disclose interests in shares | 31 |
| | UNTRACED MEMBERS | 34 |
| 72 | Power of sale | 34 |
| 73 | Application of proceeds of sale..... | 35 |
| | APPOINTMENT, RETIREMENT AND REMOVAL OF DIRECTORS | 35 |
| 74 | Number of directors | 35 |
| 75 | Power of company to appoint directors | 35 |
| 76 | Power of board to appoint directors..... | 35 |
| 77 | Appointment of executive directors | 36 |
| 78 | Eligibility for appointment as a director..... | 36 |
| 79 | Share qualification..... | 36 |
| 80 | Resolution for appointment..... | 36 |
| 81 | Re-election at intervals of no more than three years | 36 |
| 82 | Retirement of directors | 37 |
| 83 | Position of retiring director | 37 |
| 84 | Deemed re-appointment..... | 37 |
| 85 | Removal by ordinary resolution | 37 |
| 86 | Vacation of office by director..... | 38 |
| 87 | Resolution as to a vacancy conclusive..... | 39 |
| | ALTERNATE DIRECTORS | 39 |
| 88 | Appointment..... | 39 |
| 89 | Participation in board meetings..... | 39 |
| 90 | Alternate director responsible for own acts | 39 |
| 91 | Interests of alternate director | 39 |
| 92 | Revocation of appointment..... | 40 |
| | DIRECTORS' REMUNERATION, EXPENSES AND PENSIONS | 40 |

| | | |
|-----|--|----|
| 93 | Directors' fees..... | 40 |
| 94 | Expenses | 40 |
| 95 | Remuneration of executive directors..... | 41 |
| 96 | Pensions and other benefits | 41 |
| | POWERS AND DUTIES OF THE BOARD..... | 41 |
| 97 | Powers of the board | 41 |
| 98 | Powers of directors being less than minimum number | 42 |
| 99 | Powers of executive directors..... | 42 |
| 100 | Delegation to committees | 42 |
| 101 | Local management..... | 43 |
| 102 | Power of attorney | 43 |
| 103 | Use of the title "director" | 43 |
| 104 | Exercise of voting power | 44 |
| 105 | Provision for employees | 44 |
| 106 | Overseas registers | 44 |
| 107 | Borrowing powers | 44 |
| | PROCEEDINGS OF DIRECTORS AND COMMITTEES | 47 |
| 108 | Board meetings | 47 |
| 109 | Notice of board meetings..... | 47 |
| 110 | Quorum..... | 47 |
| 111 | Chairman of board..... | 48 |
| 112 | Voting..... | 48 |
| 113 | Participation by telephone or video conference..... | 48 |
| 114 | Resolution in writing | 48 |
| 115 | Proceedings of committees | 49 |
| 116 | Minutes of proceedings | 49 |
| 117 | Validity of proceedings | 50 |
| | DIRECTORS' INTERESTS..... | 50 |

| | | |
|-----|--|----|
| 118 | Director may have interests..... | 50 |
| | SEALS | 55 |
| 119 | Safe custody | 55 |
| 120 | Application of seals | 55 |
| 121 | Official seal for use abroad | 55 |
| | THE SECRETARY | 55 |
| 122 | The secretary | 55 |
| | AUTHENTICATION OF DOCUMENTS | 56 |
| 123 | Power to authenticate..... | 56 |
| | DIVIDENDS AND OTHER PAYMENTS..... | 56 |
| 124 | Declaration of dividends | 56 |
| 125 | Interim dividends | 56 |
| 126 | Entitlement to dividends | 57 |
| 127 | Calls or debts may be deducted from dividends..... | 57 |
| 128 | Distribution in specie | 57 |
| 129 | Dividends not to bear interest | 57 |
| 130 | Method of payment..... | 57 |
| 131 | Uncashed dividends | 58 |
| 132 | Unclaimed dividends | 59 |
| 133 | Payment of share dividends | 59 |
| 134 | Reserves..... | 61 |
| 135 | Capitalisation of reserves | 61 |
| 136 | Record dates..... | 63 |
| | ACCOUNTS..... | 63 |
| 137 | Inspection of records | 63 |
| 138 | Accounts to be sent to members | 63 |
| 139 | Summary financial statements..... | 64 |
| | NOTICES | 64 |

| | | |
|-----|--|----|
| 140 | Notices to be in writing | 64 |
| 141 | Service of notice on members | 64 |
| 142 | Notice in case of Death, Bankruptcy or Mental Disorder | 66 |
| 143 | Evidence of Service..... | 66 |
| 144 | Notice Binding on Transferees..... | 67 |
| 145 | Notice by Advertisement..... | 67 |
| 146 | Suspension of Postal Services | 68 |
| | WINDING UP..... | 68 |
| 147 | Division of assets | 68 |
| 148 | Transfer or sale under section 110 Insolvency Act 1986 | 68 |
| | INDEMNITY | 69 |
| 149 | Right to indemnity..... | 69 |

THE COMPANIES ACTS 1985 to 2006

PUBLIC COMPANY LIMITED BY SHARES

ARTICLES OF ASSOCIATION

OF

CYPROTEX PLC

(Amended by special resolution passed on 14 July 2009)

PRELIMINARY

1 No other regulations to apply

- 1.1 No regulations set out in any statute or in any statutory instrument or other subordinate legislation concerning companies shall apply to the company, but the following shall be the articles of association of the company.

2 Interpretation

- 2.1 In these articles, unless the context otherwise requires, the following expressions have the following meanings:

"**address**" includes any number or address used for the purposes of sending or receiving documents and/or information by electronic means;

"**articles**" means these articles of association as originally adopted or altered or varied from time to time (and "article" means one of these articles);

"**auditors**" means the auditors for the time being of the company or, in the case of joint auditors, any one of them;

"**board**" means the board of directors for the time being of the company or the directors present or deemed to be present at a duly convened meeting of directors at which a quorum is present;

"**CA 1985**" means, subject to article 2.3, the Companies Act 1985;

"**CA 2006**" means, subject to article 2.3, the Companies Act 2006;

"**chairman**" means the chairman (if any) of the board or, where the context requires, the chairman of a general meeting of the company;

"**clear days**" means (in relation to the period of a notice) that period, excluding the day when the notice is given or deemed to be given and the day for which it is given or on which it is to take effect;

"**company**" means Cyprotex plc, company number 4311107;

"**depository**" means a custodian or other person (or a nominee for such custodian or other person) appointed under contractual arrangements with the company or other arrangements approved by the board whereby such custodian or other person or nominee holds or is interested in shares of the company or rights or interests in shares of the company and issues securities or other documents of title or otherwise evidencing the entitlement of the holder thereof to or to receive such shares, rights or interests, provided and to the extent that such arrangements have been approved by the board for the purpose of these articles, and shall include, where approved by the board, the trustees (acting in their capacity as such) of any employees' share scheme established by the company or any other scheme or arrangement principally for the benefit of employees or those in the service of the company and/or its subsidiaries or their respective businesses and the managers (acting in their capacity as such) of any investment or savings plan, which in each case the board has approved;

"**director**" means a director for the time being of the company;

"**electronic copy**", "**electronic form**" and "**electronic**" have the same meanings as set out in section 1168 CA 2006;

"**execution**" includes any mode of execution (and "executed" shall be construed accordingly);

"**hard copy**" and "**hard copy form**" have the same meanings as set out in section 1168 CA 2006;

"**holder**" means (in relation to any share) the member whose name is entered in the register as the holder or, where the context permits, the members whose names are entered in the register as the joint holders, of that share;

"**London Stock Exchange**" means London Stock Exchange plc;

"**member**" means a member of the company or, where the context requires, a member of the board or of any committee;

"**office**" means the registered office for the time being of the company;

"**Operator**" has the meaning set out in the Regulations;

"**ordinary share**" means an ordinary share of £0.001 in the capital of the company;

"**paid up**" means paid up or credited as paid up;

"**recognised person**" means a recognised clearing house or a nominee of a recognised clearing house or of a recognised investment exchange which is designated as mentioned in section 778(2) CA 2006;

"**register**" means the register of members of the company to be kept pursuant to the statutes or, as the case may be, any overseas branch register kept pursuant to article 106;

"**Regulations**" means the Uncertificated Securities Regulations 2001, SI 2001 No. 3755, including any modification thereof or rules made thereunder;

"**seal**" means any common seal of the company or any securities seal that the company may be permitted to have under the statutes;

"**secretary**" means the secretary for the time being of the company or any other person appointed to perform any of the duties of the secretary of the company including (subject to the provisions of the statutes) a joint, temporary, assistant or deputy secretary;

"**securities seal**" means an official seal kept by the company pursuant to section 40 CA 1985;

"**share**" means a share in the capital of the company;

"**statutes**" means CA 1985, CA 2006 and every other statute (and any regulations subordinate thereto) for the time being in force concerning companies and affecting the company;

"**uncertificated share**" or "**participating security**" means a share which is recorded in the register as being in uncertificated form and title to which may be transferred by means of a relevant system, and references to share being held "in uncertificated form" shall be construed accordingly;

"**uncertificated proxy instruction**" means a properly authenticated dematerialised instruction (as defined in the Regulations), and/or other instruction or notification, which is sent by means of the relevant system and received by such participant in that system as the board may prescribe, in such form and subject to such terms and conditions as may from time to time be prescribed by

the board (subject always to the facilities and requirements of the relevant system concerned);

"United Kingdom" means Great Britain and Northern Ireland; and

"writing" or **"written"** means and includes printing, typewriting, lithography, photography and any other mode or modes of representing or reproducing words in a legible and non-transitory form, and includes (to the extent permitted from time to time by the statutes) electronic mail.

2.2 Unless the context otherwise requires:

- (a) words in the singular include the plural, and *vice versa*;
- (b) words importing the masculine gender include the feminine gender; and
- (c) a reference to a person includes a body corporate and an unincorporated body of persons.

2.3 A reference to any statute or provision of a statute shall include any orders, regulations or other subordinate legislation made under it and shall, unless the context otherwise requires, include any statutory modification or re-enactment of it for the time being in force.

2.4 Save as aforesaid, and unless the context otherwise requires, words or expressions contained in these articles shall bear the same meaning as in the statutes and words and expressions used in the Regulations have the same meanings when used in these articles.

2.5 The headings are inserted for convenience only and shall not affect the construction of these articles.

3 Form of resolution

3.1 Subject to the statutes, where for any purpose an ordinary resolution of the company is required, a special resolution shall also be effective.

4 Registered office

4.1 The office shall be at such place in England and Wales as the board shall from time to time appoint.

SHARE CAPITAL

5 Authorised share capital

- 5.1 The authorised share capital of the company at the date of the adoption of these articles is £300,000 divided into 300,000,000 ordinary shares of £0.001 each.

6 Uncertificated shares

- 6.1 Notwithstanding anything in these articles to the contrary, any shares in the company may be issued, held, registered, converted to, transferred or otherwise dealt with in uncertificated form and converted from uncertificated form to certificated form in accordance with the Regulations and practices instituted by the operator of the relevant system. Any provisions of these articles shall not apply to any uncertificated shares to the extent that such provisions are inconsistent with:

- (a) the holding of shares in uncertificated form;
- (b) the transfer of title to shares by means of a relevant system; or
- (c) any provision of the Regulations.

- 6.2 Without prejudice to the generality and effectiveness of the foregoing:

- (a) without prejudice to article 43 in relation to uncertificated shares, the board may also refuse to register a transfer of uncertificated shares in such other circumstances as may be permitted or required by the Regulations and the relevant system;
- (b) references in these articles to a requirement on any person to execute or deliver an instrument of transfer or certificate or other document which shall not be appropriate in the case of uncertificated shares shall, in the case of uncertificated shares, be treated as references to a requirement to comply with any relevant requirements of the relevant system and any relevant arrangements or regulations which the board may make from time to time pursuant to article 6.2(h) below;
- (c) for the purposes referred to in article 50, a person entitled by transmission to a share in uncertificated form who elects to have some other person registered shall either:
 - (i) procure that instructions are given by means of the relevant system to effect transfer of such uncertificated share to that person; or
 - (ii) change the uncertificated share to certificated form and execute an instrument of transfer of that certificated share to that person;

- (d) the company shall enter on the register the number of shares which are held by each member in uncertificated form and in certificated form and shall maintain the register in each case as is required by the Regulations and the relevant system and, unless the board otherwise determines, holdings of the same holder or joint holders in certificated form and uncertificated form shall be treated as separate holdings;
- (e) a class of share shall not be treated as two classes by virtue only of that class comprising both certificated shares and uncertificated shares or as a result of any provisions of these articles or the Regulations which applies only in respect of certificated shares or uncertificated shares;
- (f) for the purposes of article 130.1, any payment in the case of uncertificated shares may be made by means of the relevant system (subject always to the facilities and requirements of the relevant system) and without prejudice to the generality of the foregoing such payment may be made by the sending by the company or any person on its behalf of an instruction to the operator of the relevant system to credit the cash memorandum account of the holder or joint holders of such shares or, if permitted by the company, of such person as the holder or joint holders may in writing direct and for the purposes of article 130.1 the making of a payment in accordance with the facilities and requirements of the relevant system concerned shall be a good discharge to the company;
- (g) subject to the statutes the board may issue shares as certificated shares or as uncertificated shares in its absolute discretion and articles 7, 135 and 136 shall be construed accordingly;
- (h) the board may make such arrangements or regulations (if any) as it may from time to time in its absolute discretion think fit in relation to the evidencing and transfer of uncertificated shares and otherwise for the purpose of implementing and/or supplementing the provisions of this article 6 and the Regulations and the facilities and requirements of the relevant system and such arrangements and regulations (as the case may be) shall have the same effect as if set out in this article 6;
- (i) the board may utilise the relevant system to the fullest extent available from time to time in the exercise of the company's powers or functions under the statutes or these articles or otherwise in effecting any actions; and
- (j) the board may resolve that a class of shares is to become a participating security and may at any time determine that a class of shares shall cease to be a participating security.

6.3 Where any class of shares in the capital of the company is a participating security and the company is entitled under any provisions of the statutes or the rules made and practices instituted by the Operator of any relevant system or under these articles to dispose of, forfeit, enforce a lien or sell or otherwise procure the sale of any shares which are held in uncertificated form, such entitlement (to the extent permitted by the Regulations and the rules made and practices instituted by the Operator of the relevant system) shall include the right to:

- (a) request or require the deletion of any computer-based entries in the relevant system relating to the holding of such shares in uncertificated form; and/or
- (b) require any holder of any uncertificated shares which are the subject of any exercise by the company of any such entitlement, by notice in writing to the holder concerned, to change his holding of such uncertificated shares into certificated form within such period as may be specified in the notice, prior to completion of any disposal, sale or transfer of such shares or direct the holder to take such steps, by instructions given by means of a relevant system or otherwise, as may be necessary to sell or transfer such shares; and/or
- (c) send a notification in writing to the Operator requiring conversion of those shares into certificated form (such conversion being required to enable the company to deal with such shares in accordance with the articles); and/or
- (d) appoint or require the holder of any shares to appoint by notice any person to take such other steps, by instruction given by means of a relevant system or otherwise, in the name of the holder of such shares as may be required to effect transfer of such shares and such steps shall be as effective as if they had been taken by the registered holder of the uncertificated shares concerned; and/or
- (e) transfer any uncertificated shares which are the subject of any exercise by the company of any such entitlement by entering the name of the transferee in the register in respect of that share as a transferred share; and/or
- (f) otherwise rectify or change the register in respect of that share in such manner as may be appropriate; and
- (g) take such other action as may be necessary to enable those shares to be registered in the name of the person to whom the shares have been sold or disposed of or as directed by him.

6.4 For the purposes of this article 6:

- (a) words and expressions shall have the same respective meanings as in the Regulations;
- (b) references herein to an uncertificated share or to a share (or to a holding of shares) being in uncertificated form are references to that share being an uncertificated unit of a security, and references to a certificated share or to a share being in certificated form are references to that share being a unit of a security which is not an uncertificated unit; and
- (c) "cash memorandum account" means an account so designated by the Operator of the relevant system.

7 Allotment

- 7.1 Subject to the provisions of the statutes and to any relevant authority of the company in general meeting required by the statutes all unissued shares created in the capital of the company shall be at the disposal of the board, which may allot (with or without conferring rights of renunciation), grant options over, offer or otherwise deal with or dispose of them or rights to subscribe for or convert any security into shares to such persons (including the directors themselves), at such times and generally on such terms and conditions as the board may decide, provided that no share shall be issued at a discount.

8 Redeemable shares

- 8.1 Subject to the provisions of the statutes and to any special rights for the time being attached to any existing shares, any share may be issued which is, or at the option of the company or of the holder of such share is liable, to be redeemed on such terms and in such manner as these articles may provide.

9 Power to attach rights

- 9.1 Subject to the provisions of the statutes and to any special rights for the time being attached to any existing shares, any shares may be allotted or issued with, or have attached to them, such preferred, deferred or other special rights or restrictions, whether in regard to dividend, voting, transfer, return of capital or otherwise, as the company may from time to time by ordinary resolution determine or, if no such resolution has been passed or so far as the resolution does not make specific provision, as the board may determine.

10 Share warrants to bearer

- 10.1 The company may, with respect to any fully paid shares, issue a warrant ("**share warrant**") stating that the bearer of the warrant is entitled to the shares specified in it and may provide (by coupons or otherwise) for the payment of future dividends on the shares included in a share warrant.

10.2 The powers referred to in article 10.1 may be exercised by the board, which may determine and vary the conditions on which share warrants shall be issued and, in particular, on which:

- (a) a new share warrant or coupon will be issued in the place of one damaged, defaced, worn out or lost (provided that no new share warrant shall be issued to replace one that has been lost unless the board is satisfied beyond reasonable doubt that the original has been destroyed);
- (b) the bearer of a share warrant shall be entitled to receive notice of and to attend, vote and demand a poll at general meetings;
- (c) dividends will be paid; and
- (d) a share warrant may be surrendered and the name of the holder entered in the register in respect of the shares specified in it.

Subject to such conditions and to these articles, the bearer of a share warrant shall be deemed to be a member for all purposes. The bearer of a share warrant shall be subject to the conditions for the time being in force and applicable thereto, whether made before or after the issue of such share warrant.

11 Commission and brokerage

11.1 The company may, in connection with the issue of any shares, exercise all powers of paying commission and brokerage conferred or permitted by the statutes. Subject to the provisions of the statutes, any such commission or brokerage may be satisfied by the payment of cash, the allotment of fully or partly paid shares, the grant of an option to call for an allotment of shares or any combination of such methods.

12 Trusts not to be recognised

12.1 Except as otherwise expressly provided by these articles, as required by law or as ordered by a court of competent jurisdiction, the company shall not recognise any person as holding any share on any trust, and (except as aforesaid) the company shall not be bound by or recognise (even if having notice of it) any equitable, contingent, future, partial or other claim to or interest in any share except an absolute right of the holder to the whole of the share.

ALTERATION OF SHARE CAPITAL

13 Increase, consolidation, cancellation and sub-division

13.1 The company may from time to time by ordinary resolution:

- (a) increase its share capital by such sum to be divided into shares of such amount as the resolution prescribes;

- (b) consolidate and divide all or any of its share capital into shares of a larger nominal amount than its existing shares;
- (c) cancel any shares which at the date of the passing of the resolution have not been taken or agreed to be taken by any person, and reduce the amount of its share capital by the amount of the shares so cancelled; and
- (d) subject to the provisions of the statutes, sub-divide its shares or any of them into shares of smaller nominal amount, and may by such resolution determine that, as between the shares resulting from such sub-division, one or more of the shares may, as compared with the others, have any such preferred, deferred or other special rights or be subject to any such restrictions as the company has power to attach to unissued or new shares.

14 Fractions of shares

14.1 Whenever, as the result of any consolidation, division or sub-division of shares, any difficulty arises, the board may settle it as it thinks fit, and, in particular (but without prejudice to the generality of the foregoing), where the number of shares held by any holder is not an exact multiple of the number of shares to be consolidated into a single share and as a result of such consolidation such holder would become entitled to a fraction of a consolidated share:

- (a) the board may determine which of the shares of such holder are to be treated as giving rise to such fractional entitlement and may decide that any of those shares shall be consolidated with any of the shares of any other holder or holders which are similarly determined by it to be treated as giving rise to a fractional entitlement for such other holder or holders, into a single consolidated share and the board may, on behalf of all such holders, sell such consolidated share for the best price reasonably obtained to any person (including the company) and distribute the net proceeds of sale after deduction of the expenses of sale in due proportion among those holders (except that any amount otherwise due to a holder, being less than £1.00 or such other sum as the board may from time to time determine, may be retained for the benefit of the company); or
- (b) provided that the necessary unissued shares are available, the board may issue to such holder credited as fully paid by way of capitalisation the minimum number of shares required to round up his holding to an exact multiple of the number of shares to be consolidated into a single share (such issue being deemed to have been effected prior to consolidation); and the amount required to pay up such shares shall be appropriated at the board's discretion from any of the sums standing to the credit of any of the company's reserve accounts (including share premium account and capital redemption reserve) or to the credit of profit and loss account and

capitalised by applying the same in paying up the share. In relation to such a capitalisation, the board may exercise all the powers conferred on it by article 135 without an ordinary resolution of the company.

- 14.2 For the purposes of any sale of consolidated shares pursuant to article 14.1, the board may authorise some person to execute an instrument of transfer (or, as the case may be, to procure a transfer pursuant to the rules of a relevant system) of the shares to, or in accordance with the directions of, the purchaser, and the transferee shall not be bound to see to the application of the purchase money, nor shall his title to the shares be affected by any irregularity in or invalidity of the proceedings in reference to the sale.

15 Reduction of capital

- 15.1 Subject to the provisions of the statutes, these articles and to any rights for the time being attached to any shares, the company may by special resolution reduce its share capital or any capital redemption reserve or share premium account or other undistributable reserve in any way.

16 Purchase of own shares

- 16.1 Subject to the provisions of the statutes, these articles and any rights for the time being attached to any shares, the company may purchase any of its own shares of any class (including any redeemable shares). Any shares to be so purchased may be selected in any manner whatsoever.

VARIATION OF CLASS RIGHTS

17 Sanction to variation

- 17.1 If at any time the share capital of the company is divided into shares of different classes, the special rights attached to any class may be varied or abrogated by a special resolution passed at a separate meeting of the holders of the relevant class of shares in accordance with CA 2006.

18 Class meetings

- 18.1 All the provisions in these articles as to general meetings shall mutatis mutandis apply to every meeting of the holders of any class of shares. The board may convene a meeting of the holders of any class of shares whenever it thinks fit and whether or not the business to be transacted involves a variation or abrogation of class rights. The quorum at every such meeting shall be not less than two persons holding or representing by proxy at least one-third of the nominal amount paid up on the issued shares of the class. Every holder of shares of the class, present in person or by proxy, may demand a poll. Each such holder shall on a poll be entitled to one vote for every share of the class held by him. If at any adjourned meeting of such holders such quorum as

aforesaid is not present, not less than one person holding shares of the class who is present in person or by proxy shall be a quorum.

19 Deemed variation

- 19.1 Subject to the terms of issue of or rights attached to any shares, the rights or privileges attached to any class of shares shall be deemed not to be varied or abrogated by the creation or issue of any new shares ranking *pari passu* in all respects (save as to the date from which such new shares shall rank for dividend) with or subsequent to those already issued or by the reduction of the capital paid up on such shares or by the purchase or redemption by the company of its own shares in accordance with the provisions of the statutes and these articles.

EVIDENCE OF TITLE

20 Right to share certificates

- 20.1 On becoming the holder of any certificated share, every person (except a recognised person in respect of whom the company is not by law required to complete and have ready for delivery a certificate) shall be entitled, without charge, to have issued within two months after allotment or lodgement of a transfer (unless the terms of issue of the shares provide otherwise) one certificate for all the certificated shares of each class registered in his name. Such certificate shall specify the number, class and distinguishing numbers (if any) of the shares in respect of which it is issued and the amount or respective amounts paid up thereon and shall be issued as provided in article 120.
- 20.2 The issued shares of a particular class which are fully paid up and rank *pari passu* for all purposes shall not bear a distinguishing number. All other shares shall bear a distinguishing number.
- 20.3 The company shall not be bound to issue more than one certificate in respect of shares held jointly by two or more persons. Delivery of a certificate to the person first named on the register as the holder of such shares shall be sufficient delivery to all joint holders.
- 20.4 Where a member (other than a recognised person) has transferred part only of the shares comprised in a certificate, he shall be entitled to a certificate for the balance of such shares without charge.
- 20.5 No certificate shall be issued representing shares of more than one class or in respect of shares held by a recognised person, or in respect of any uncertificated shares.

21 Replacement certificates

- 21.1 Any two or more certificates representing certificated shares of any one class held by any member may, at his request, be cancelled and a single new certificate for such shares issued in lieu without charge on surrender of the original certificates for cancellation.
- 21.2 If any member shall surrender for cancellation a share certificate representing certificated shares held by him and request the company to issue in lieu two or more share certificates representing such shares in such proportions as he may specify, the board may, if it thinks fit, comply with such request.
- 21.3 Share certificates may be renewed or replaced on such terms as to provision of evidence and indemnity (with or without security) and to payment of any exceptional out of pocket expenses, including those incurred by the company in investigating such evidence and preparing such indemnity and security, as the board may decide, and on surrender of the original certificate (where it is defaced, damaged or worn out), but without any further charge.
- 21.4 In the case of shares held jointly by several persons, any such request as is mentioned in this article 21 may be made by any one of the joint holders.

LIEN ON SHARES

22 Lien on shares not fully paid

- 22.1 The company shall have a first and paramount lien on each of its shares which is not fully paid, for all amounts payable to the company (whether presently or not) in respect of that share and to the extent and in the circumstances permitted by the statutes. The company's lien on such a share shall extend to every amount (including, without limitation, dividends). The board may waive any lien which has arisen and may resolve that any share shall for some limited period be exempt wholly or partially from the provisions of this article.

23 Enforcement of lien by sale

- 23.1 The board may sell all or any of the shares subject to any lien at such time or times and in such manner as it may determine. However, no sale shall be made until such time as the moneys in respect of which such lien exists or some part thereof are or is presently payable or the liability or engagement in respect of which such lien exists is liable to be presently fulfilled or discharged, and until a demand and notice in writing stating the amount due or specifying the liability or engagement and demanding payment or fulfilment or discharge thereof and giving notice of intention to sell in default shall have been served on the holder or the persons (if any) entitled by transmission to the shares, and default in payment, fulfilment or discharge shall have been made by him or them for fourteen clear days after service of such notice. For giving effect to

any such sale, the board may authorise some person to execute an instrument of transfer (or, as the case may be, to procure a transfer pursuant to the rules of a relevant system) of the shares sold in the name and on behalf of the holder or the persons entitled by transmission in favour of the purchaser or as the purchaser may direct. The purchaser shall not be bound to see to the application of the purchase money, and the title of the transferee to the shares shall not be affected by any irregularity in or invalidity of the proceedings in reference to the sale.

24 Application of proceeds of sale

- 24.1 The net proceeds of any sale of shares subject to any lien, after payment of the costs, shall be applied in or towards satisfaction of so much of the amount due to the company or of the liability or engagement (as the case may be) as is presently payable or is liable to be presently fulfilled or discharged. The balance (if any) shall (in the case of a sale of certificated shares, on surrender to the company for cancellation of the certificate for the shares sold), subject to a like lien for any moneys not presently payable or any liability or engagement not liable to be presently fulfilled or discharged as existed on the shares before the sale, be paid to the holder or the person (if any) entitled by transmission to the shares so sold, without interest.

CALLS ON SHARES

25 Calls

- 25.1 Subject to the terms of allotment of shares, the board may from time to time make calls on the members in respect of any moneys unpaid on the shares, of any class, held by them respectively (whether in respect of nominal value or premium) and not payable on a date fixed by or in accordance with the terms of issue. Each member shall (subject to receiving at least fourteen clear days' written notice specifying when and where payment is to be made and whether or not by instalments) be liable to pay the amount of every call so made on him as required by the notice. A call shall be deemed to have been made at the time when the resolution of the board authorising such call was passed or (as the case may require) any person to whom power has been delegated pursuant to these articles serves notice of exercise of such power. A call may be required to be paid by instalments and may, before receipt by the company of any sum due thereunder, be either revoked or postponed in whole or part as regards all or any such members as the board may determine. A person on whom a call is made shall remain liable notwithstanding the subsequent transfer of the shares in respect of which the call was made.

26 Liability of joint holders

- 26.1 The joint holders of a share shall be jointly and severally liable for the payment of all calls in respect thereof.

27 Interest on calls

- 27.1 If the whole of the sum payable in respect of any call is not paid on or before the day appointed for payment, the person from whom it is due and payable shall pay all costs, charges and expenses that the company may have incurred by reason of such non-payment, together with interest on the unpaid amount from the day appointed for payment thereof to the time of actual payment at the rate fixed by the terms of the allotment of the share or in the notice of the call or, if no rate is so fixed, at such rate, not exceeding ten per cent per annum (compounded on a six monthly basis), as the board shall determine. The board may waive payment of such costs, charges, expenses or interest in whole or in part.

28 Rights of member when call unpaid

- 28.1 Unless the board otherwise determines, no member shall be entitled to receive any dividend or to be present or vote or speak at a general meeting or at any separate general meeting of the holders of any class of shares either in person or (save as proxy for another member) by proxy, or be reckoned in a quorum, or to exercise any other right or privilege as a member in respect of a share held by him unless and until he shall have paid all calls for the time being due and payable by him in respect of that share, whether alone or jointly with any other person, together with interest and expenses (if any) to the company.

29 Sums due on allotment treated as calls

- 29.1 Any sum payable in respect of a share on allotment or at any fixed date, whether in respect of the nominal value of the share or by way of premium or as an instalment of a call, shall, for all purposes of these articles be deemed to be a call duly made. If it is not paid, the provisions of these articles shall apply as if such amount had become due and payable by virtue of a call.

30 Power to differentiate

- 30.1 The board may make arrangements on the allotment or issue of shares for a difference as between the allottees or holders of such shares in the amount and time of payment of calls.

31 Payment in advance of calls

- 31.1 The board may, if it thinks fit, receive from any member willing to advance the same all or any part of the moneys uncalled and unpaid on the shares held by him. Such payment in advance of calls shall extinguish *pro tanto* the liability on

the shares on which it is made. The company may pay interest on the money paid in advance, or so much of it as exceeds the amount for the time being called up on the shares in respect of which such advance has been made, at such rate as the board may decide. The board may at any time repay the amount so advanced on giving to such member not less than three months' notice in writing of its intention in that behalf, unless before the expiration of such notice the amount so advanced shall have been called up on the shares in respect of which it was advanced.

32 Delegation of power to make calls

- 32.1 If any uncalled capital of the company is included in or charged by any mortgage or other security, the board may delegate on such terms as it thinks fit to the person in whose favour such mortgage or security is executed, or to any other person in trust for him, the power to make calls on the members in respect of such uncalled capital, to sue in the name of the company or otherwise for the recovery of moneys becoming due in respect of calls so made and to give valid receipts for such moneys. The power so delegated shall subsist during the continuance of the mortgage or security, notwithstanding any change of directors, and shall be assignable if expressed so to be.

FORFEITURE OF SHARES

33 Notice if call not paid

- 33.1 If any member fails to pay the whole of any call or any instalment of any call on or before the day appointed for payment, the board may at any time serve a notice in writing on such member or on any person entitled to the shares by transmission, requiring payment, on a date not less than fourteen clear days from the date of the notice, of the amount unpaid and any interest which may have accrued thereon and any costs, charges and expenses incurred by the company by reason of such non-payment. The notice shall name the place where the payment is to be made and state that, if the notice is not complied with, the shares in respect of which such call was made will be liable to be forfeited.

34 Forfeiture for non-compliance

- 34.1 If the notice referred to in article 33 is not complied with, any share in respect of which it was given may, at any time before the payment required by the notice has been made, be forfeited by a resolution of the board to that effect. Such forfeiture shall include all dividends declared or other moneys payable in respect of the forfeited shares and not paid before the forfeiture.

35 Notice after forfeiture

- 35.1 When any share has been forfeited, notice of the forfeiture shall be served on the person who was before forfeiture the holder of the share or the person entitled to

such share by transmission (as the case may be). An entry of such notice having been given, and of the forfeiture with the date thereof, shall forthwith be made in the register in respect of such share. However, no forfeiture shall be invalidated by any omission to give such notice or to make such entry as aforesaid.

36 Forfeiture may be annulled

36.1 The board may, at any time before any share so forfeited has been cancelled or sold, re-allotted or otherwise disposed of, annul the forfeiture, on terms that payment shall be made of all calls and interest due thereon and all expenses incurred in respect of the share and on such further terms (if any) as the board shall see fit.

37 Surrender

37.1 The board may accept a surrender of any share liable to be forfeited. In such case, references in these articles to forfeiture shall include surrender.

38 Disposal of forfeited shares

38.1 Every share which shall be forfeited shall thereupon become the property of the company. Subject to the provisions of the statutes, any such share may be sold, re-allotted or otherwise disposed of, either to the person who was before forfeiture the holder thereof or entitled thereto or to any other person, on such terms and in such manner as the board shall determine. The board may, for the purposes of the disposal, authorise some person to transfer the share in question and may enter the name of the transferee in respect of the transferred share in the register notwithstanding (in the case of a certificated share) the absence of any share certificate being lodged in respect thereof. An instrument of transfer executed by that person, or a transfer effected by that person pursuant to the rules of a relevant system, shall be as effective as if it had been executed or effected by the holder of, or the person entitled by transmission to, the share. The company may receive the consideration (if any) given for the share on its disposal.

39 Effect of forfeiture

39.1 A member whose shares have been forfeited shall cease to be a member in respect of the shares forfeited and shall surrender to the company for cancellation the certificate (if any) for such shares. He shall nevertheless be liable to pay to the company all calls made and not paid on such shares at the time of forfeiture, and interest thereon from the date of the forfeiture to the date of payment, in the same manner in all respects as if the shares had not been forfeited, and to satisfy all (if any) claims, demands and liabilities which the company might have enforced in respect of the shares at the time of forfeiture, without any reduction or allowance for the value of the shares at the time of forfeiture or for any consideration received on their disposal.

40 Extinction of claims

40.1 The forfeiture of a share shall involve the extinction at the time of forfeiture of all interest in and all claims and demands against the company in respect of the share and all other rights and liabilities incidental to the share as between the holder whose share is forfeited and the company, except only such of those rights and liabilities as are by these articles expressly saved, or as are by the statutes given or imposed in the case of past members.

41 Evidence of forfeiture

41.1 A statutory declaration by a director or the secretary that a share has been forfeited in pursuance of these articles, and stating the date on which it was forfeited, shall, as against all persons claiming to be entitled to the share adversely to the forfeiture thereof, be conclusive evidence of the facts therein stated. The declaration, together with the receipt of the company for the consideration (if any) given for the share on the sale or disposition thereof shall (subject, if necessary, to the execution of an instrument of transfer or the transfer of the share pursuant to the rules of a relevant system) constitute a good title to the share. Subject to the completion of any necessary transfer, such person shall be registered as the holder of the share and shall be discharged from all calls made prior to such sale or disposition and shall not be bound to see to the application of the purchase money or other consideration (if any), nor shall his title to the share be affected by any act, omission or irregularity relating to or connected with the proceedings in reference to the forfeiture or disposal of the share. Such person shall not (except by express agreement with the company) become entitled to any dividend which might have accrued on the share before the completion of the sale or disposition thereof.

TRANSFER OF SHARES

42 Transfer of certificated shares

42.1 Subject to such of the restrictions of these articles as may be applicable, each member may transfer all or any of his certificated shares by instrument of transfer in writing in any usual form or in any form approved by the board. Such instrument shall be executed by or on behalf of the transferor and (in the case of a transfer of a certificated share which is not fully paid up) by or on behalf of the transferee. The transferor shall be deemed to remain the holder of such share until the name of the transferee is entered in the register in respect of it. All instruments of transfer which are registered may be retained by the company.

43 Transfer of uncertificated shares

43.1 Subject to such of the restrictions of these articles as may be applicable, each member may transfer all or any of his uncertificated shares in accordance with the

rules and regulations in force from time to time which are applicable to a relevant system.

44 Right to refuse registration

44.1 The board may in its absolute discretion and without giving any reason, refuse to register any transfer of a share (or renunciation of a renounceable letter of allotment) unless:

- (a) it is in respect of a share which is fully paid up;
- (b) it is in respect of only one class of shares;
- (c) it is in favour of a single transferee or not more than four joint transferees;
- (d) it is duly stamped (if so required);
- (e) (if it is in respect of a certificated share) it is delivered for registration to the office or such other place as the board may from time to time determine, accompanied (except in the case of a transfer by a recognised person where a certificate has not been issued or in the case of a renunciation) by the certificate for the shares to which it relates and such other evidence as the board may reasonably require to prove the title of the transferor or person renouncing and the due execution of the transfer or renunciation by him or, if the transfer or renunciation is executed by some other person on his behalf, the authority of that person to do so; and
- (f) in the case of an uncertificated share, in the circumstances as set out in the Regulations.

Notwithstanding the provisions of this article 44, provided that the board shall not refuse to register any transfer or renunciation of partly paid shares which are admitted to trading on the Alternative Investment Market of the London Stock Exchange on the grounds that they are partly paid shares in circumstances where such refusal would prevent dealings in such shares from taking place on an open and proper basis.

45 Notice of refusal

If the board refuses to register a transfer of a share, it shall, within two months after the date on which the transfer was lodged with the company, send notice of the refusal to the transferee together with the reason(s) for the refusal. Any instrument of transfer which the board refuses to register shall (except in the case of suspected or actual fraud) be returned to the person depositing it.

46 Closing of register

46.1 The registration of transfers of shares or of any class of shares may be suspended at such times and for such periods (not exceeding thirty days in any year) as the board may from time to time determine, subject to the board having first obtained any consent or authorisation to such closure which may be required. Notice of closure of the register shall be given in accordance with the requirements of the statutes.

47 Fees on registration

47.1 No fee shall be charged for registration of a transfer or on the registration of any probate, letters of administration, certificate of death or marriage, power of attorney, notice or other instrument relating to or affecting the title to any shares.

48 Transfers by renunciation

48.1 Nothing in these articles shall preclude the board from recognising a renunciation of the allotment of any share by the allottee in favour of some other person.

TRANSMISSION OF SHARES

49 On death

49.1 If a member dies, the survivors or survivor, where he was a joint holder, and his executors or administrators, where he was a sole or the only survivor of joint holders, shall be the only persons recognised by the company as having any title to his shares. Nothing in these articles shall release the estate of a deceased member from any liability in respect of any share which has been solely or jointly held by him.

50 Election of person entitled by transmission

50.1 Any person becoming entitled to a share in consequence of the death or bankruptcy of any member, or of any other event giving rise to a transmission of such entitlement by operation of law, may, on such evidence as to his title being produced as the board may require, elect either to become registered as a member or to have some person nominated by him registered as a member. If he elects to become registered himself, he shall give notice to the company to that effect. If he elects to have some other person registered, he shall execute an instrument of transfer of such share to that person or, as the case may be, procure the transfer of such share to that person pursuant to the rules of a relevant system. All the provisions of these articles relating to the transfer of shares shall apply to the notice or transfer (as the case may be) as if it were effected by the member and his death, bankruptcy or other event as aforesaid had not occurred. Where the entitlement of a person to a share in consequence of the death or bankruptcy of a member or of any other event giving rise to its transmission by

operation of law is proved to the satisfaction of the board, the board shall within two months after proof cause the entitlement of that person to be noted in the register.

51 Rights on transmission

- 51.1 Where a person becomes entitled to a share in consequence of the death or bankruptcy of any member, or of any other event giving rise to a transmission of such entitlement by operation of law, the rights of the holder in relation to such share shall cease. However, the person so entitled may give a good discharge for any dividends and other moneys payable in respect of it and shall have the same rights to which he would be entitled if he were the holder of the share, except that he shall not, before he is registered as the holder of the share, be entitled in respect of it to receive notice of, or to attend or vote at, any meeting of the company or at any separate meeting of the holders of any class of shares of the company. The board may at any time give notice requiring any such person to elect either to be registered himself or to transfer the share. If the notice is not complied with within sixty days, the board may thereafter withhold payment of all dividends and other moneys payable in respect of such share until the requirements of the notice have been complied with.

DESTRUCTION OF DOCUMENTS

52 Destruction of documents

- 52.1 The company may destroy:
- (a) any instrument of transfer, after six years from the date on which it is registered;
 - (b) any dividend mandate or any variation or cancellation thereof or any notification of change of name or address, after two years from the date on which it is recorded;
 - (c) any share certificate, after one year from the date on which it is cancelled; and
 - (d) any other document on the basis of which any entry in the register is made, after six years from the date on which an entry was first made in the register in respect of it,

provided that the company may destroy any such type of document at a date earlier than that authorised by this article if a copy of such document is retained on microfilm or by other similar means on which such copy is retained until the expiration of the period applicable to the destruction of the original of such document.

- 52.2 It shall be conclusively presumed in favour of the company that every entry in the register purporting to have been made on the basis of a document so destroyed was duly and properly made, that every instrument of transfer so destroyed was duly registered, that every share certificate so destroyed was duly cancelled, that every other document so destroyed had been properly dealt with in accordance with its terms and was valid and effective in accordance with the particulars in the records of the company, provided that:
- (a) this article 52 shall apply only to the destruction of a document in good faith and without notice of any claim (regardless of the parties to it) to which the document might be relevant;
 - (b) nothing in this article 52 shall be construed as imposing on the company any liability in respect of the destruction of any such document otherwise than as provided for in this article 52 which would not attach to the company in the absence of this article 52; and
 - (c) references in this article 52 to the destruction of any document include references to the disposal of it in any manner.

GENERAL MEETINGS

53 Annual general meetings

- 53.1 Subject to the provisions of the statutes, annual general meetings shall be held at such time and place as the board may determine.

54 Convening of general meetings

- 54.1 The board may convene a general meeting whenever it thinks fit. A general meeting shall also be convened on such requisition, or in default may be convened by such requisitionists, as provided by section 303 CA 2006. At any meeting convened on such requisition or by such requisitionists no business shall be transacted except that stated by the requisition or proposed by the board. If there are not within the United Kingdom sufficient members of the board to convene a general meeting, any director may call a general meeting.

55 Notice of general meetings

- 55.1 A notice of a general meeting shall specify:
- (a) whether the meeting is an annual general meeting or a general meeting;
 - (b) the place, the day and the time of the meeting;
 - (c) the general nature of the business of the meeting;

- (d) if the meeting is convened to consider a special resolution, the intention to propose the resolution as such; and
 - (e) with reasonable prominence, that a member entitled to attend and vote is entitled to appoint one or more proxies to attend, speak and vote instead of him and that a proxy need not also be a member.
- 55.2 The notice shall be given to the members (other than any who, under the provisions of these articles or of any restrictions imposed on any shares, are not entitled to receive notice from the company), to the directors and to the auditors.
- 55.3 If, after a notice convening a general meeting of the company has been despatched to members, the board becomes aware of any fact, event or circumstances which, in the board's opinion, would make it impractical or inappropriate to hold the general meeting on the date or time or at the place for which notice has been given, the board may give notice pursuant to article 145 to those entitled to receive the notice pursuant to article 55.2 either cancelling such meeting or postponing such meeting to a time and date which is not less than twenty one clear days from the date of such notice, which shall also specify the place at which such postponed meeting shall be held.
- 55.4 The board may specify in the notice of meeting a time by which a person must be entered on the register in order to have the right to attend or vote at the meeting. The time specified must not be more than 48 hours before the time fixed for the meeting.
- 55.5 If, after sending a notice of meeting but before the meeting is held, or after the adjournment of a meeting but before the adjourned meeting is held (whether or not notice of the adjourned meeting is required), the board decides that it is impractical or unreasonable, for a reason beyond its control, to hold the meeting at the declared place and/or time, it may change the place and/or postpone the time at which the meeting is held. If such a decision is made, the board may then change the place and/or postpone the time again if it decides that it is reasonable to do so. In either case:
- (a) no new notice of the meeting need be sent, but the board shall, if practicable, advertise the new place and/or date and time of the meeting in at least two newspapers having a national circulation and shall make arrangements for notice of the change of place and/or postponement of date and time to appear at the original place and/or date and time; and
 - (b) a proxy appointment in relation to the meeting may be received at the address specified in the notice of meeting not less than 48 hours before any postponed time appointed for holding the meeting.

56 Omission to send notice

- 56.1 The accidental omission to send a notice of meeting or, in cases where it is intended that it be sent out with the notice, an instrument of proxy to, or the non-receipt of either by, any person entitled to receive the same shall not invalidate the proceedings at that meeting.

PROCEEDINGS AT GENERAL MEETINGS

57 Quorum

- 57.1 No business shall be transacted at any annual general meeting or general meeting unless a quorum is present when the meeting proceeds to business. Two qualifying persons entitled to vote on the business to be transacted shall be a quorum unless :

- (a) each is a qualifying person only because he is authorised to act as the representative of a corporation in relation to the meeting and they are representatives of the same corporation; or
- (b) each is a qualifying person only because he is appointed as a proxy of a member in relation to the meeting and they are proxies of the same member.

- 57.2 For the purposes of this article 57, a "qualifying person" means:

- (a) an individual who is a member of the company;
- (b) a person authorised to act as the representative in relation to the meeting of a corporation which is a member of the company; or
- (c) a person appointed in relation to the meeting as a proxy of a member of the company.

58 If quorum not present

- 58.1 If within thirty minutes (or such longer interval as the chairman in his absolute discretion thinks fit) from the time appointed for the holding of an annual general meeting or a general meeting a quorum is not present, or if during a meeting such a quorum ceases to be present, the meeting, if convened on the requisition of members, shall be dissolved. In any other case, the meeting shall stand adjourned to such time (not being less than seven nor more than sixty days later) and place as the chairman (or, in default, the board) may determine. If at such adjourned meeting a quorum is not present within fifteen minutes from the time appointed for holding the meeting, the meeting shall be dissolved.

59 Chairman

- 59.1 The chairman of the board shall preside at every annual general meeting and general meeting of the company. If there be no such chairman, or if at any meeting he shall not be present within five minutes after the time appointed for holding the meeting, or shall be unwilling to act as chairman, the deputy chairman (if any), if present and willing to do so shall preside at such meeting, but if neither the chairman or deputy chairman is present and willing to act, the directors present shall choose one of their number to act as chairman or, if there be only one director present, he shall be chairman if willing to act. If there be no director present and willing to act, the members present and entitled to vote shall choose one of their number to be chairman of the meeting.
- 59.2 The chairman of the meeting who presides pursuant to the provisions of article 59.1 may, at any time during a general meeting of the company, nominate any director of the company to be the chairman of the meeting for the remainder of or for any part of the meeting.

60 Directors may attend and speak

- 60.1 A director shall, notwithstanding that he is not a member, be entitled to attend and speak at any annual general meeting or general meeting and at any separate meeting of the holders of any class of shares of the company.

61 Power to adjourn

- 61.1 The chairman may, with the consent of a meeting at which a quorum is present, and shall, if so directed by the meeting, adjourn any meeting from time to time (or indefinitely) and from place to place as the meeting shall determine. However, without prejudice to any other power which he may have under these articles or at common law, the chairman may, without the need for the consent of the meeting, interrupt or adjourn any meeting from time to time and from place to place or for an indefinite period if he is of the opinion that it has become necessary to do so in order to secure the comfort, safety and security of those attending and the proper and orderly conduct of the meeting or to give all persons entitled to do so a reasonable opportunity of attending, speaking and voting at the meeting or to ensure that the business of the meeting is properly disposed of.

62 Notice of adjourned meeting

- 62.1 Where a meeting is adjourned indefinitely, the board shall fix the time and place for the adjourned meeting. Whenever a meeting is adjourned for fourteen days or more or indefinitely, seven clear days' notice at the least, specifying the place, the day and time of the adjourned meeting and the general nature of the business to be transacted, shall be given in the same manner as in the case of an original meeting. Save as aforesaid, no member shall be entitled to any notice of an adjournment or of the business to be transacted at any adjourned meeting.

63 Business of adjourned meeting

63.1 No business shall be transacted at any adjourned meeting other than the business which might properly have been transacted at the meeting from which the adjournment took place.

64 Accommodation of members and security arrangements

64.1 The board may, for the purpose of controlling the level of attendance and ensuring the safety and security of those attending at any place specified for the holding of any annual general meeting or general meeting, from time to time make such arrangements as the board shall consider to be appropriate in the circumstances and may from time to time vary any such arrangements or make new arrangements in place thereof. In the case of any meeting to which such arrangements apply the board may, when specifying the place of the meeting:

- (a) direct that the meeting shall be held at a place specified in the notice at which the chairman of the meeting shall preside ("**the principal place**"); and
- (b) make arrangements for simultaneous attendance and participation at other places by members otherwise entitled to attend the general meeting but who cannot be accommodated in the principal place and who are excluded therefrom under the provisions of this article or who wish to attend at any of such other places, provided that persons attending at the principal place and at any of such other places shall be able to see, and hear and be seen and heard by, persons attending at the principal place and at such other places, by any means.

Such arrangements for simultaneous attendance may include arrangements for regulating the level of attendance in any manner aforesaid as between the principal place and any of such other places, provided that they shall operate so that any member who cannot be accommodated in the principal place as aforesaid are able to attend at one of such other places. For the purposes of all other provisions of these articles any such meeting shall be treated as being held and taking place at the principal place.

64.2 For the purpose of ensuring the safety and security of those attending any meeting the board may require that any person wishing to attend any meeting should submit to such searches or other security arrangements as the board shall consider appropriate in the circumstances and shall be entitled in its absolute discretion to refuse entry to any meeting to any person who refuses to cooperate with or to submit to such searches or to otherwise comply with such security arrangements.

VOTING

65 Method of voting

65.1 At any general meeting, a resolution put to a vote of the meeting shall be decided on a show of hands, unless (before or on the declaration of the result of the show of hands) a poll is duly demanded. Subject to the provisions of the statutes, a poll may be demanded by:

- (a) the chairman of the meeting; or
- (b) by at least five members present in person or by proxy and entitled to vote at the meeting; or
- (c) a member or members present in person or by proxy representing not less than one-tenth of the total voting rights of all the members having the right to vote at the meeting; or
- (d) a member or members present in person or by proxy holding shares conferring a right to vote at the meeting, being shares on which an aggregate sum has been paid up equal to not less than one-tenth of the total sum paid up on all the shares conferring that right.

65.2 The chairman of the meeting may also demand a poll before a resolution is put to the vote on a show of hands;

65.3 At annual general meetings and general meetings, resolutions shall be put to the vote by the chairman and there shall be no requirement for the resolution to be proposed or seconded by any person.

66 Chairman's declaration conclusive on show of hands

66.1 Unless a poll is duly demanded and the demand is not withdrawn, a declaration by the chairman of the meeting that a resolution has on a show of hands been carried, or carried unanimously or by a particular majority, or lost, or not carried by a particular majority, and an entry to that effect in the book containing the minutes of proceedings of the company, shall be conclusive evidence thereof, without proof of the number or proportion of the votes recorded in favour of or against such resolution.

67 Objection to error in voting

67.1 No objection shall be raised to the qualification of any voter or to the counting of, or failure to count, any vote, except at the meeting or adjourned meeting at which the vote objected to is given or tendered or at which the error occurs. Any objection or error shall be referred to the chairman of the meeting and shall only vitiate the decision of the meeting on any resolution if the chairman

decides that the same is of sufficient magnitude to vitiate the resolution or may otherwise have affected the decision of the meeting. The decision of the chairman on such matters shall be final and conclusive.

68 Amendment to resolutions

- 68.1 If an amendment shall be proposed to any resolution under consideration but shall in good faith be ruled out of order by the chairman of the meeting, any error in such ruling shall not invalidate the proceedings on the substantive resolution.
- 68.2 In the case of a resolution duly proposed as a special resolution, no amendment thereto (other than an amendment to correct a patent error) may in any event be considered or voted on and, in the case of a resolution duly proposed as an ordinary resolution, no amendment thereto (other than an amendment to correct a patent error) may be considered or voted on unless either, at least forty-eight hours prior to the time appointed for holding the meeting or adjourned meeting at which such ordinary resolution is to be proposed, notice in writing of the terms of the amendment and intention to move the same has been lodged at the office or received in electronic form at such address (if any) for the time being specified by or on behalf of the company for that purpose (or the address that the company has been deemed to agree by the statutes) or the chairman of the meeting in his absolute discretion decides that it may be considered or voted on.
- 68.3 With the consent of the chairman of the meeting, an amendment to a resolution may be withdrawn by its proposer before that resolution is put to the vote.

69 Procedure on a poll

- 69.1 Any poll duly demanded on the election of a chairman of a meeting or on any question of adjournment shall be taken forthwith. A poll duly demanded on any other matter shall be taken at such time and place, not being more than thirty days from the date of the meeting or adjourned meeting at which the poll was demanded, as the chairman shall direct. The chairman shall determine the manner (including the use of ballot or voting papers or tickets or electronic mail) in which a poll shall be taken and may appoint scrutineers who need not be members. No notice need be given of a poll not taken immediately if the time and place at which it is to be taken are announced at the meeting at which it is demanded. In any other case, at least seven clear days' notice shall be given specifying the time and place at which the poll is to be taken. The result of the poll shall be deemed to be the resolution of the meeting at which the poll was demanded.
- 69.2 The demand for a poll (other than on the election of a chairman or any question of adjournment) shall not prevent the continuance of the meeting for the transaction of any business other than the question on which a poll has been demanded. If a

poll is demanded before the declaration of the result on a show of hands and the demand is duly withdrawn, the meeting shall continue as if the demand had not been made.

- 69.3 The demand for a poll may, before the poll is taken, be withdrawn, but only with the consent of the chairman. A demand for a poll so withdrawn shall validate the result of a show of hands declared before the demand was made.
- 69.4 On a poll, votes may be given in person or by proxy. A member entitled to more than one vote need not, if he votes, use all his votes or cast all the votes he uses in the same way.

70 Votes of members

- 70.1 Subject to any special rights or restrictions as to voting attached to any shares by or in accordance with these articles, on a show of hands every member who is present in person or by proxy or (being a corporation) present by a duly authorised representative shall have one vote and on a poll every member who is present in person or by proxy or (being a corporation) present by a duly authorised representative shall have one vote for every share of which he is the holder.
- 70.2 In the case of joint holders of a share, the vote of the senior who tenders a vote, whether in person or by proxy, shall be accepted to the exclusion of the votes of the other joint holders, and for this purpose seniority shall be determined by the order in which the names of the holders stand in the register in respect of the share.
- 70.3 No member shall, unless the directors otherwise determine, be entitled in respect of shares held by him to vote at any annual general meeting or general meeting or at any separate meeting of the holders of any class of shares in the company either personally or by proxy, or to exercise any privilege as a member, unless all calls or other sums presently payable by him in respect of shares in the company have been paid.
- 70.4 Where in England or elsewhere a receiver or other person (by whatever name called) has been appointed by any court claiming jurisdiction in that regard to exercise powers with respect to the property or affairs of any member on the ground (however formulated) of mental disorder, the directors may in their absolute discretion, upon or subject to production of such evidence of the appointment as the directors may require, permit such receiver or other person on behalf of such member to vote in person or by proxy at any general meeting or to exercise any other right conferred by membership in relation to meetings of the company.
- 70.5 No objection shall be raised to the admissibility of any vote except at the meeting or adjourned meeting at which the vote objected to is or may be given

or tendered, and every vote not disallowed at such meeting shall be valid for all purposes. Any such objection made in due time shall be referred to the chairman of the meeting whose decision shall be final and conclusive.

70.6 On a poll votes may be given either personally or by proxy. On a poll a member entitled to more than one vote need not, if he votes, use all his votes or cast all the votes he uses in the same way.

70.7 Any person (whether a member or not) may be appointed to act as a proxy. The appointment of a proxy (whether made by an instrument in writing or by electronic means), shall be in any usual or common form, or such other form as may be approved by the directors, and shall be signed by the appointor or by his attorney duly authorised in writing, or if the appointor is a corporation shall be either under its common seal or under the hand of a duly authorised officer or attorney of the corporation. The directors may, but shall not be bound to, require evidence of the authority of such officer or attorney. The appointment of a proxy need not be witnessed.

70.8 The appointment of a proxy together with (unless the directors waive such requirement) the power of attorney or other authority (if any) under which it is signed, or a notarially certified or office copy of such power or authority, shall in the case of an instrument in writing be deposited, at the office, or at such other place in the United Kingdom as may be specified for that purpose in the notice calling the meeting or in any instrument of proxy sent out by the company in relation to the meeting, not less than 48 hours before the time appointed for holding the meeting or adjourned meeting at which the person named in the instrument proposes to vote, and in default the instrument of proxy shall not be treated as valid. The like time limit shall also apply to the cancellation or revocation of any such instrument. In the case of an appointment of a proxy by electronic means, where an address has been specified for the purpose of receiving communications by electronic means in the notice convening the meeting, or in any instrument of proxy sent out by the company in relation to the meeting, or in any invitation contained in a communication in electronic form to appoint a proxy issued by the company in relation to the meeting, such appointment must be received at such address not less than 48 hours before the time for holding the meeting or adjourned meeting at which the person named in the appointment proposes to vote. Subject to article 70.2, where more than one appointment of a proxy is delivered, deposited or received in respect of the same shares, that delivered, deposited or received latest shall prevail; if it is not clear which was delivered, deposited or received latest, none shall be valid. The appointment of a proxy to vote at any meeting delivered, deposited or received as previously stated shall, unless the contrary is stated on it, be valid to empower the proxy so appointed to vote on any poll taken or demanded at such meeting or at any adjournment of such meeting. The appointment of a proxy shall not be valid after the expiration of 12 months from the date named in it as the date of its execution, except on a poll demanded at

a meeting or an adjourned meeting in cases where the original meeting was held within 12 months from such date. In calculating the periods mentioned in this article, the directors may specify, in any case, that no account shall be taken of any part of a day that is not a working day.

- 70.9 The appointment of a proxy shall be deemed to include the right to demand or join in demanding a poll but shall not confer any further right to speak at the meeting, except with the permission of the chairman of the meeting.
- 70.10 A vote given in accordance with the terms of the appointment of a proxy or by the duly authorised representative of a corporate member, or a poll demanded by proxy or by the duly authorised representative of a corporate member, shall be valid notwithstanding (in the case of a proxy) the previous death or mental disorder of the principal or the revocation of the appointment of a proxy or of the authority under which the appointment of a proxy was executed or (in the case of a duly authorised representative of a corporate member) the revocation of his appointment, provided that no intimation in writing or electronic form of such death, mental disorder or revocation shall have been received by the company at the office (or, where the appointment of the proxy was contained in a communication by electronic means, at the address at which such appointment was duly received) or at such other place at which it was required to be delivered, deposited or received under article 70.8 at least three hours before the commencement of the meeting or adjourned meeting at which the vote is given or the poll demanded or (in the case of a poll taken otherwise than on the same day as the meeting or adjourned meeting) the time appointed for taking the poll.

71 Failure to disclose interests in shares

- 71.1 If a member, or any other person appearing to be interested in shares held by that member, has been issued with a notice pursuant to section 793 CA 2006 (a "**section 793 notice**") and has failed in relation to any shares ("**the default shares**", which expression includes any shares issued after the date of such notice in respect of those shares) to give the company the information thereby required within the prescribed period from the service of the notice, the following sanctions shall apply unless the board otherwise determines:
- (a) the member shall not be entitled in respect of the default shares to be present or to vote (either in person or by representative or proxy) at any annual general meeting or general meeting or at any separate meeting of the holders of any class of shares or on any poll or to exercise any other right conferred by membership in relation to any such meeting or poll; and
 - (b) where the default shares represent at least 0.25 per cent in nominal value of the issued shares of their class:

- (i) any dividend or other money payable in respect of the shares shall be withheld by the company, which shall not have any obligation to pay interest on it, and the member shall not be entitled to elect, pursuant to article 133, to receive shares instead of that dividend; and
- (ii) (subject to the requirements of the relevant system where the default shares are held in uncertificated form) no transfer, other than an excepted transfer, of any default shares held by the member shall be registered unless:
 - (A) the member is not himself in default as regards supplying the information required; and
 - (B) the member proves to the satisfaction of the board that no person in default as regards supplying such information is interested in any of the shares the subject of the transfer.

71.2 Where the sanctions under article 71.2 apply in relation to any default shares, they shall cease to have effect (and any dividends withheld under article 71.2(b) shall become payable):

- (a) if the shares are transferred by means of an excepted transfer, but only in respect of the shares transferred; or
- (b) at the end of the period of seven days (or such shorter period as the board may determine) following receipt by the company of the information required by the notice mentioned in that paragraph and the board being fully satisfied that such information is full and complete.

71.3 Where, on the basis of information obtained from a member in respect of any share held by him, the company issues a section 793 notice to any other person, it shall at the same time send a copy of the notice to the member, but the accidental omission to do so, or the non-receipt by the member of the copy, shall not invalidate or otherwise affect the application of article 71.1.

71.4 Where default shares in which a person appears to be interested are held by a depositary, the provisions of this article 71 shall be treated as applying only to those shares held by the depositary in which such person appears to be interested and not (insofar as such person's apparent interest is concerned) to any other shares held by the depositary.

71.5 Where the member on which a section 793 notice is served is a depositary acting in its capacity as such, the obligations of the depositary as a member of the company shall be limited to disclosing to the company such information relating to any person appearing to be interested in the shares held by it as has been

recorded by it pursuant to the arrangements entered into by the company or approved by the board pursuant to which it was appointed as a depositary.

71.6 For the purposes of this article 71:

- (a) a person, other than the member holding a share, shall be treated as appearing to be interested in that share if the member has informed the company that the person is, or may be, so interested or if the company (after taking account of any information obtained from the member or, pursuant to a notice under section 793 CA 2006, from anyone else) knows or has reasonable cause to believe that the person is, or may be, so interested;
- (b) "interested" shall be construed as it is for the purpose of section 821 CA 2006;
- (c) reference to a person having failed to give the company the information required by a section 793 notice, or being in default as regards supplying such information, includes reference:
 - (i) to his having failed or refused to give all or any part of it; and
 - (ii) to his having given information which he knows to be false in a material particular, or having recklessly given information which is false in a material particular;
- (d) "prescribed period" means fourteen days from the date of service from the section 793 notice;
- (e) "excepted transfer" means, in relation to any shares held by a member:
 - (i) a transfer by way of, or pursuant to, acceptance of a takeover offer for the company (within the meaning of section 974 CA 2006); or
 - (ii) a transfer in consequence of a sale made through a recognised investment exchange (as defined in section 285 of the Financial Services and Markets Act 2000) or any other stock exchange outside the United Kingdom on which the company's shares are normally traded; or
 - (iii) a transfer which is shown to the satisfaction of the board to be made in consequence of a sale of the whole of the beneficial interest in the shares to a person who is unconnected with the member and with any other person appearing to be interested in the shares.

71.7 Nothing contained in this article 71 shall be taken to limit the powers of the company under section 794 CA 2006.

UNTRACED MEMBERS

72 Power of sale

72.1 The company shall be entitled to sell at the best price reasonably obtainable any share of a member, or any share to which a person is entitled by transmission, if and provided that:

- (a) during the period of twelve years prior to the date of the publication of the advertisements referred to in sub-paragraph (b) below (or, if published on different dates, the earlier or earliest thereof) no cheque, order or warrant in respect of such share sent by the company through the post in a pre-paid envelope addressed to the member or to the person entitled by transmission to the share, at his address on the register or other last known address given by the member or person to which cheques, orders or warrants in respect of such share are to be sent has been cashed and the company has received no communications in respect of such share from such member or person, provided that during such period of twelve years the company has paid at least three cash dividends (whether interim or final) and no such dividend has been claimed by the person entitled to it;
- (b) on or after expiry of the said period of twelve years the company has given notice of its intention to sell such share by advertisements in two newspapers of which one shall be a national newspaper published in the United Kingdom and the other shall be a newspaper circulating in the area of the address on the register or other last known address of the member or the person entitled by transmission to the share or the address for the service of notices notified under article 141.4;
- (c) the said advertisements, if not published on the same day, shall have been published within thirty days of each other; and
- (d) during the further period of three months following the date of publication of the said advertisements (or, if published on different dates, the later or latest thereof) and prior to the exercise of the power of sale the company has not received any communication in respect of such share from the member or person entitled by transmission.

72.2 To give effect to any sale of shares pursuant to this article, the board may authorise some person to transfer the shares in question and may enter the name of the transferee in respect of the transferred shares in the register notwithstanding the absence of any share certificate being lodged in respect thereof and may issue a new certificate to the transferee. An instrument of

transfer executed (or, as the case may be, a transfer effected pursuant to the rules of a relevant system) by that person shall be as effective as if it had been executed or effected by the holder of, or the person entitled by transmission to, the shares. The purchaser shall not be bound to see to the application of the purchase moneys, nor shall his title to the shares be affected by any irregularity or invalidity in the proceedings relating to the sale.

- 72.3 If during the period of twelve years referred to in article 72.1, or during any period ending on the date when all the requirements of paragraphs (a) to (d) of article 72.1 have been satisfied, any additional shares have been issued in respect of those held at the beginning of, or previously so issued during, any such period and all the requirements of paragraphs (b) to (d) of article 72.1 have been satisfied in regard to such additional shares, the company shall also be entitled to sell the additional shares.

73 Application of proceeds of sale

- 73.1 The company shall account to the member or other person entitled to such share for the net proceeds of such sale by carrying all moneys in respect thereof to a separate account. The company shall be deemed to be a debtor to, and not a trustee for, such member or other person in respect of such moneys. Moneys carried to such separate account may either be employed in the business of the company or invested in such investments as the board may from time to time think fit. No interest shall be payable to such member or other person in respect of such moneys and the company shall not be required to account for any money earned on them.

APPOINTMENT, RETIREMENT AND REMOVAL OF DIRECTORS

74 Number of directors

- 74.1 The number of directors shall not be less than two or more than fifteen.

75 Power of company to appoint directors

- 75.1 Subject to the provisions of these articles, the company may by ordinary resolution appoint a person who is willing to act to be a director, either to fill a vacancy or as an addition to the existing board.

76 Power of board to appoint directors

- 76.1 Without prejudice to the power of the company to appoint any person to be a director pursuant to these articles, the board shall have power at any time to appoint any person who is willing to act as a director, either to fill a vacancy or as an addition to the existing board. Any director so appointed shall retire and be eligible for election at the annual general meeting of the company next following

such appointment and shall not be taken into account in determining the number of directors who are to retire by rotation at such meeting.

77 Appointment of executive directors

77.1 Subject to the provisions of the statutes, the board may from time to time appoint one or more of its body to hold any employment or executive office (including that of executive chairman, chief executive or managing director) for such term (subject to the provisions of the statutes) and subject to such other conditions as the board thinks fit in accordance with article 99. The board may revoke or terminate any such appointment without prejudice to any claim for damages for breach of contract between the director and the company.

78 Eligibility for appointment as a director

78.1 No person, other than a director retiring (by rotation or otherwise), shall be appointed or re-appointed a director at any annual general meeting or general meeting unless:

- (a) he is recommended by the board; or
- (b) not less than seven nor more than forty-two clear days before the date appointed for the meeting, notice duly executed by a member (other than the person to be proposed) entitled to vote at the meeting has been given to the company of the intention to propose that person for appointment or re-appointment, stating the particulars which would, if he were so appointed or re-appointed, be required to be included in the company's register of directors, together with notice executed by that person of his willingness to be appointed or re-appointed, is lodged at the office.

79 Share qualification

79.1 A director shall not be required to hold any shares of the company.

80 Resolution for appointment

80.1 A resolution for the appointment of two or more persons as directors by a single resolution shall be void unless an ordinary resolution that it shall be so proposed has first been agreed to by the meeting without any vote being given against it.

81 Re-election at intervals of no more than three years

81.1 A director will retire from office at the third annual general meeting following his appointment or following his last re-appointment by members at an annual general meeting.

82 Retirement of directors

82.1 Subject to article 82.2, at each annual general meeting of the company, all those directors who have been in office for three years or more since their election or last re-election shall retire from office. In addition, any director who wishes to retire from office may do so at any annual general meeting.

82.2 At each annual general meeting of the company, all those non-executive directors who have been in office for nine years or more since the date on which they were originally elected as a non-executive director shall retire from office.

83 Position of retiring director

83.1 A director who retires at an annual general meeting (whether by rotation or otherwise) may, if willing to act, be re-appointed. If he is not re-appointed or deemed to have been re-appointed, he shall retain office until the meeting appoints someone in his place or, if it does not do so, until the end of the meeting.

84 Deemed re-appointment

84.1 At any annual general meeting at which a director retires under any provision of these articles the company may by ordinary resolution fill the office being vacated by electing thereto the retiring director or some other person eligible for election. In default, the retiring director shall be deemed to have been elected except in any of the following cases:

- (a) where at such meeting it is expressly resolved not to fill such office or a resolution for the election of such director is put to the meeting and lost;
- (b) where such director has given notice in writing to the company that he is unwilling to be re-elected;
- (c) where the default is due to the moving of a resolution in contravention of article 80; or
- (d) where such director is for any reason no longer eligible for re-election.

85 Removal by ordinary resolution

85.1 In addition to any power of removal conferred by the statutes, the company may by ordinary resolution remove any director before the expiration of his period of office, but without prejudice to any claim for damages which he may have for breach of any contract of service between him and the company, and may (subject to these articles) by ordinary resolution appoint another person who is willing to act to be a director in his place. Any person so appointed shall be treated, for the purposes of determining the time at which he or any other director

is to retire, as if he had become a director on the day on which the person in whose place he is appointed was last appointed or re-appointed a director.

86 Vacation of office by director

86.1 Without prejudice to the provisions for retirement of directors contained in these articles, the office of a director shall be vacated if:

- (a) he resigns by notice in writing delivered to the secretary at the office or tendered at a board meeting;
- (b) he ceases to be a director by virtue of any provision of the statutes, is removed from office pursuant to these articles or the statutes or becomes prohibited by law from being a director;
- (c) he becomes bankrupt, has an interim receiving order made against him, makes any arrangement or compounds with his creditors generally or applies to the court for an interim order under section 253 of the Insolvency Act 1986 in connection with a voluntary arrangement under that Act;
- (d) an order is made by any court of competent jurisdiction on the ground (howsoever formulated) of mental disorder for his detention or for the appointment of a guardian or receiver or other person to exercise powers with respect to his affairs or he is admitted to hospital in pursuance of an application for admission for treatment under the Mental Health Act 1983 (as amended) or, in Scotland, under the Mental Health (Care and Treatment) (Scotland) Act 2003 and the board resolves that his office be vacated;
- (e) both he and his alternate director appointed pursuant to the provisions of these articles (if any) are absent, without the permission of the board, from board meetings for six consecutive months and the board resolves that his office be vacated;
- (f) he is removed from office by notice in writing signed by all of the other directors (without prejudice to any claim for damages which he may have for breach of any contract between him and the company) and, for this purpose, a set of like notices each signed by, one or more of the directors shall be as effective as a single notice signed by the requisite number of directors; or
- (g) in the case of any director who holds any executive office with the company, his appointment as such is terminated or expires and the directors resolve that his office be vacated.

87 Resolution as to a vacancy conclusive

- 87.1 A resolution of the board declaring a director to have vacated office under the terms of article 86 shall be conclusive as to the fact and grounds of vacation stated in the resolution.

ALTERNATE DIRECTORS

88 Appointment

- 88.1 Each director (other than an alternate director) may, by notice in writing to the company in hard copy form or electronic form sent or delivered to such address (if any) specified by the company for that purpose or, in default of such specification, to the office, appoint any other director or any person approved for that purpose by the board and willing to act, to be his alternate.
- 88.2 No appointment of an alternate director who is not already a director shall be effective until his consent to act as a director in the form prescribed by the statutes has been received at the office.
- 88.3 An alternate director need not hold a share qualification and shall not be counted in reckoning any maximum or minimum number of directors allowed by these articles.

89 Participation in board meetings

- 89.1 Every alternate director shall (subject to his giving to the company an address within the United Kingdom at which notices may be served on him) be entitled to receive notice of all meetings of the board and all committees of the board of which his appointor is a member and, in the absence from such meetings of his appointor, to attend and vote at such meetings and to exercise all the powers, rights, duties and authorities of his appointor. A director acting as alternate director shall have a separate vote at board meetings for each director for whom he acts as alternate director, but he shall count as only one for the purpose of determining whether a quorum is present.

90 Alternate director responsible for own acts

- 90.1 Every person acting as an alternate director shall be an officer of the company, shall alone be responsible to the company for his own acts and defaults and shall not be deemed to be the agent of the director appointing him.

91 Interests of alternate director

- 91.1 An alternate director shall save as otherwise provided in these articles be entitled to contract and be interested in and benefit from contracts or arrangements with the company and to be repaid expenses and to be indemnified to the same extent

mutatis mutandis as if he were a director. However, he shall not be entitled to receive from the company any fees for his services as alternate, except only such part (if any) of the fee payable to his appointor as such appointor may by notice in writing to the company direct. Subject to this article, the company shall pay to an alternate director such expenses as might properly have been paid to him if he had been a director and to indemnify him to the same extent as if he were a director.

92 Revocation of appointment

92.1 An alternate director shall cease to be an alternate director:

- (a) if his appointor revokes his appointment; or
- (b) if his appointor ceases for any reason to be a director, provided that if any director retires but is re-appointed or deemed to be re-appointed at the same meeting, any valid appointment of an alternate director which was in force immediately before his retirement shall remain in force; or
- (c) if any event happens in relation to him which, if he were a director otherwise appointed, would cause him to vacate office.

DIRECTORS' REMUNERATION, EXPENSES AND PENSIONS

93 Directors' fees

93.1 The maximum aggregate annual fees payable to the directors (excluding any director holding salaried employment in the company or any subsidiary of the company) for services in the office of director shall be the sum of £250,000 or such larger sum as the company by ordinary resolution shall from time to time determine. Any fees payable pursuant to this article shall be distinct from any salary, remuneration or other amounts payable to a director pursuant to any other provisions of these articles and shall accrue from day to day.

94 Expenses

94.1 Each director shall be entitled to be repaid all reasonable travelling, hotel and other expenses properly incurred by him in or about the performance of his duties as a director, including any expenses incurred in attending meetings of the board or any committee of the board or general meetings or separate meetings of the holders of any class of shares or of debentures of the company. If in the opinion of the directors it is desirable that any of their number should go or reside abroad or make any special journeys or perform any special services on behalf of the company or its business, such director or directors may be paid such reasonable additional remuneration by way of salary, percentage of profits or otherwise and expenses therefor as the directors may from time to time determine.

95 Remuneration of executive directors

95.1 The salary or remuneration of any director appointed to hold any employment or executive office in accordance with the provisions of these articles may be either a fixed sum of money, or may altogether or in part be governed by business done or profits made or otherwise determined by the board, and may be in addition to or in lieu of any fee payable to him for his services as director pursuant to these articles.

96 Pensions and other benefits

96.1 The board may exercise all the powers of the company to provide pensions or other retirement or superannuation benefits and to provide death or disability benefits or other allowances or gratuities (whether by insurance or otherwise) for, or to institute and maintain any institution, association, society, club, trust, other establishment or profit-sharing, share incentive, share purchase or employees' share scheme calculated to advance the interests of the company or to benefit, any person who is or has at any time been a director or employee of the company or any company which is a holding company or a subsidiary undertaking of or allied to or associated with the company or any such holding company or subsidiary undertaking or any predecessor in business of the company or of any such holding company or subsidiary undertaking, and for any member of his family (including a spouse or former spouse) and any person who is or was dependent on him. For such purpose the board may establish, maintain, subscribe and contribute to any scheme, institution, association, club, trust or fund and pay premiums and, subject to the provisions of the statutes, lend money or make payments to, guarantee or give an indemnity in respect of, or give any financial or other assistance in connection with any of the aforesaid matters. The board may procure any of such matters to be done by the company either alone or in conjunction with any other person. Any director or former director shall be entitled to receive and retain for his own benefit any pension or other benefit provided under this article and shall not be obliged to account for it to the company.

POWERS AND DUTIES OF THE BOARD

97 Powers of the board

97.1 Subject to the provisions of the statutes, the memorandum of association of the company and these articles and to any directions given by special resolution of the company, the business of the company shall be managed by the board, which may exercise all the powers of the company, whether relating to the management of the business or not. No alteration of the memorandum of association or of these articles and no such direction given by the company shall invalidate any prior act of the board which would have been valid if such alteration had not been made or such direction had not been given. Provisions contained elsewhere in these

articles as to any specific power of the board shall not be deemed to limit the general powers given by this article.

98 Powers of directors being less than minimum number

98.1 If the number of directors is less than the minimum for the time being prescribed by these articles, the remaining director or directors shall act only for the purposes of appointing an additional director or directors to make up such minimum or of convening a general meeting of the company for the purpose of making such appointment. If no director or directors are able or willing to act, any two members may summon a general meeting for the purpose of appointing directors. Subject to the provisions of these articles, any additional director so appointed shall hold office only until the dissolution of the annual general meeting of the company next following such appointment unless he is re-elected during such meeting.

99 Powers of executive directors

99.1 The board may from time to time:

- (a) delegate or entrust to and confer on any director holding executive office (including a chief executive or managing director) such of its powers, authorities and discretions (with power to sub-delegate) for such time, on such terms and subject to such conditions as it thinks fit; and
- (b) revoke, withdraw, alter or vary all or any of such powers.

100 Delegation to committees

100.1 The board may delegate any of its powers, authorities and discretions (with power to sub-delegate) for such time on such terms and subject to such conditions as it thinks fit to any committee consisting of one or more directors and (if thought fit) one or more other persons, provided that:

- (a) a majority of the members of a committee shall be directors; and
- (b) no resolution of a committee shall be effective unless a majority of those present when it is passed are directors or alternate directors.

100.2 The board may confer such powers either collaterally with, or to the exclusion of and in substitution for, all or any of the powers of the board in that respect and may from time to time revoke, withdraw, alter or vary any of such powers and discharge any such committee in whole or in part. Insofar as any power, authority or discretion is so delegated, any reference in these articles to the exercise by the board of such power, authority or discretion shall be construed as if it were a reference to the exercise of such power, authority or discretion by such committee.

101 Local management

101.1 The board may establish any local or divisional boards or agencies for managing any of the affairs of the company in any specified locality, either in the United Kingdom or elsewhere, and may appoint any persons to be members of such local or divisional board, or any managers or agents, and may fix their remuneration. The board may delegate to any local or divisional board, manager or agent so appointed any of its powers, authorities and discretions (with power to sub-delegate) and may authorise the members for the time being of any such local or divisional board, or any of them, to fill any vacancies and to act notwithstanding vacancies; and any such appointment or delegation may be made for such time, on such terms and subject to such conditions as the board may think fit. The board may confer such powers either collaterally with, or to the exclusion of and in substitution for, all or any of the powers of the board in that respect and may from time to time revoke, withdraw, alter or vary all or any of such powers. Subject to any terms and conditions expressly imposed by the board, the proceedings of any local or divisional board or agency with two or more members shall be governed by such of these articles as regulate the proceedings of the board, so far as they are capable of applying.

102 Power of attorney

102.1 The board may, by power of attorney or otherwise, appoint any person or persons to be the agent of the company and may delegate to any such person or persons any of its powers, authorities and discretions (with power to sub-delegate), in each case for such purposes and for such time, on such terms (including as to remuneration) and subject to such conditions as it thinks fit (including any provisions for the protection and convenience of any person dealing with the agent). The board may confer such powers either collaterally with, or to the exclusion of and in substitution for, all or any of the powers of the board in that respect and may from time to time revoke, withdraw, alter or vary any of such powers.

103 Use of the title "director"

103.1 The board may appoint any person (not being a director) to any office or employment having a designation or title including the word "director" or attach to any existing office or employment with the company such designation or title and may terminate any such appointment or the use of such designation or title. The inclusion of the word "director" in the designation or title of any such office or employment shall not imply that such person is, or is deemed to be, or is empowered in any respect to act as, a director for any of the purposes of the statutes or these articles.

104 Exercise of voting power

104.1 The board may exercise or cause to be exercised the voting power conferred by the shares in any other company held or owned by the company, or any power of appointment to be exercised by the company, in such manner in all respects as it thinks fit (including the exercise of the voting power or power of appointment in favour of the appointment of any director as a director or other officer or employee of such company or in favour of the payment of remuneration to the directors, officers or employees of such company).

105 Provision for employees

105.1 Pursuant to section 719 CA 1985, to the extent that it is in force from time to time, and section 247 CA 2006 from the time that it comes into force, the directors are hereby authorised to make such provision as may seem appropriate for the benefit of any persons employed or formerly employed by the company or any of its subsidiary undertakings in connection with the cessation or the transfer of the whole or part of the undertaking of the company or any subsidiary undertaking. Any such provision shall be made by a resolution of the directors in accordance with such sections.

106 Overseas registers

106.1 Subject to the provisions of the statutes, the board may exercise the powers conferred on the company with regard to the keeping of an overseas branch, local or other register and may make and vary such regulations as it thinks fit respecting the keeping of any such register.

107 Borrowing powers

107.1 The board may exercise all the powers of the company to borrow money and to mortgage or charge all or any part of the undertaking, property and assets (present and future) and uncalled capital of the company and, subject to the provisions of the statutes, to create and issue debenture and other loan stock and debentures and other securities, whether outright or as collateral security for any debt, liability or obligation of the company or of any third party.

107.2 The board shall restrict the borrowings of the company and exercise all voting and other rights and powers of control exercisable by the company in respect of its subsidiary undertakings so as to procure (as regards its subsidiary undertakings in so far as it can procure by such exercise) that the aggregate principal amount at any one time outstanding in respect of moneys borrowed by the group (exclusive of moneys borrowed by one group company from another and after deducting cash deposited) shall not, at any time, without the previous sanction of an ordinary resolution of the company, exceed an amount equal to:

- (i) the nominal amount of the share capital of the company issued and paid up or credited as paid up; and
- (ii) the consolidated reserves of the group.

107.3 For the purpose of this article 107:

- (a) the expression "the group" means the company and its subsidiaries;
- (b) the expression "moneys borrowed" includes the following except in so far as otherwise taken into account under this article:
 - (i) the nominal or principal amount of any share capital debentures or borrowed moneys (together in each case with any fixed or minimum premium payable on the final date for repayment) of any person the beneficial interest wherein is not for the time being owned by a member of the group and the repayment whereof is guaranteed by a member of the group;
 - (ii) the principal amount (together with any fixed or minimum premium payable on the final date for repayment) of any debenture (whether secured or unsecured) of a member of the group owned otherwise than by another member of the group;
 - (iii) the principal amount raised by acceptances under any acceptance credit granted by any bank or accepting house;
- (c) there shall not be taken into account as "moneys borrowed":
 - (i) monies borrowed by a member of the group for the purpose of repaying or otherwise discharging other moneys brought into account and so to be applied within four months of being so borrowed pending their application for such purpose within such period;
 - (ii) that proportion of the total monies borrowed of any partly-owned subsidiary of the company which its issued equity share capital not for the time being beneficially owned directly or indirectly by the company bears to the whole of its issued equity share capital but a like proportion of any borrowings from such subsidiary by the company shall fall to be treated as borrowings of the company concerned notwithstanding the same would not otherwise fall to be taken into account; and
 - (iii) borrowings of any company becoming a subsidiary of the company for a period of six months from the date of its becoming a

subsidiary to the extent that such borrowings do not exceed its borrowings outstanding on the date of its becoming a subsidiary;

- (d) there shall be offset against the amount of any monies borrowed by the company and its subsidiaries, subject, in the case of any held or deposited by a partly-owned subsidiary, to the exclusion therefrom of a proportion equal to the proportion of the equity share capital of the partly-owned subsidiary which is not attributable directly or indirectly to the company, an amount equal to the aggregate of all cash in hand, credit balances on current or deposit account with banks, cash deposits, certificates of deposit and debt securities of governments and companies and similar instruments owned by the company and/or any of its subsidiaries which are or represent amounts available for repayment of any monies borrowed falling to be taken into account for the purposes of sub-paragraph (b)(i) – (iii) above of this article;
- (e) the expression "the consolidated reserves of the group" means the amounts standing to the credit of the reserves of the group including any share premium account, capital redemption reserve fund, taxation equalisation account and profit and loss account, but excluding:
 - (i) any amounts attributable to goodwill and other intangible assets;
 - (ii) the amount of any debit balance of profit and loss account;
 - (iii) any amount set aside for taxation other than tax equalisation;
 - (iv) amounts attributable to minority members;
- (f) the paid-up share capital of the company and the consolidated reserves of the group at any time shall be taken to be:
 - (i) those shown in the latest audited consolidated balance sheet then available but adjusted as may be necessary in respect of (1) changes in the paid-up share capital of the company since the date of that balance sheet and (2) any company which since the date of such balance sheet has become or has ceased to be a subsidiary of the company; or
 - (ii) pending an audited consolidated balance sheet being available, the shares of the company in issue on incorporation but adjusted as may be necessary in respect of (1) changes in the paid-up share capital of the company since the date of its incorporation and (2) any company which since the date of the company's incorporation has become or has ceased to be a subsidiary of the company.

107.4 No debt incurred or security given in respect of moneys borrowed in excess of the limit imposed by this article 107 shall be invalid or ineffectual, except in the case of express notice to the lender or recipient of the security at the time when the debt was incurred or security given that the limit had been or would thereby be exceeded, but no lender or other person dealing with the company shall be concerned to see or enquire whether such limit is observed.

PROCEEDINGS OF DIRECTORS AND COMMITTEES

108 Board meetings

108.1 Subject to the provisions of these articles, the board may meet for the despatch of business, adjourn and otherwise regulate its proceedings as it thinks fit.

109 Notice of board meetings

109.1 One director may, and the secretary at the request of a director shall, summon a board meeting at any time on reasonable notice. Notice of a board meeting shall be deemed to be properly given to a director if it is given to him personally or by word of mouth or (subject to article 109.2) sent in writing to him in hard copy form or electronic form at such address as he may from time to time specify for this purpose or if he does not specify such address at his last known address or any other address given by him to the company for that purpose. A director may waive the requirement that notice be given to him of any board meeting, either prospectively or retrospectively.

109.2 When summoning a board meeting, it shall be necessary to use reasonable endeavours to give notice of that meeting to a director who is absent from the United Kingdom.

109.3 A director who is absent from the United Kingdom shall not, in any event, be entitled to any longer period of notice than if he had been present in the United Kingdom. The accidental failure to give notice to, or the non-receipt of notice by, any director of any board meeting shall not invalidate the proceedings at that board meeting.

110 Quorum

110.1 The quorum necessary for the transaction of business may be determined by the board and until otherwise determined shall be three directors save where the company has two directors or one director only in which case the quorum shall be two directors or one director as the case may be. A duly convened meeting of the board at which a quorum is present shall be competent to exercise all or any of the authorities, powers, and discretions for the time being vested in or exercisable by the board.

111 Chairman of board

111.1 The board may appoint one of its body chairman and one of its body deputy chairman of its meetings and may determine the period for which they are to hold office and may at any time remove either or both of them from office. If no such chairman or deputy chairman is elected, or if at any meeting neither a chairman nor a deputy chairman is present within five minutes of the time appointed for holding the same, the directors present shall choose one of their number to be chairman of such meeting. Any chairman or deputy chairman may also hold executive office under the company.

112 Voting

112.1 Questions arising at any board meeting shall be determined by a majority of votes. In the case of an equality of votes, the chairman of that meeting shall have a second or casting vote.

113 Participation by telephone or video conference

113.1 Any director or his alternate may validly participate in a meeting of the board or a committee of the board through the medium of conference telephone, video conferencing link or any other form of electronic means, provided that all persons participating in the meeting are able to hear and speak to each other throughout such meeting, or by a series of telephone calls or video conferences or by exchange of communications in any other electronic form from and to the chairman of the meeting.

113.2 A person so participating by such means as set out in article 113.1 with the chairman of the meeting shall be deemed to be present in person at the meeting and shall accordingly be counted in a quorum and be entitled to vote. Such a meeting shall be deemed to take place where the largest group of those participating is assembled or, if there is no group which is larger than any other group, where the chairman of the meeting then is.

113.3 A resolution passed at any meeting held in the above manner, and signed by the chairman of the meeting, shall be as valid and effectual as if it had been passed at a meeting of the board (or committee, as the case may be) duly convened and held.

114 Resolution in writing

114.1 A resolution in writing signed by all the directors for the time being entitled to receive notice of a board meeting and to vote on the resolution at a board meeting and not being less than a quorum, or by all the members of a committee of the board for the time being entitled to receive notice of such committee meeting and to vote on the resolution at such committee meeting and not being less than a quorum of that committee, shall be as valid and effective for all purposes as a

resolution duly passed at a meeting of the board (or committee, as the case may be). For this purpose:

- (a) a director signifies his agreement to a proposed written resolution when the company receives from him a document indicating his agreement to the resolution authenticated in the manner permitted by the statutes for a document in the relevant form;
- (b) a director so signifying his agreement to a proposed written resolution may send the document in hard copy form or in electronic form to such address (if any) for the time being specified by the company for that purpose or, in default of such specification, to the office;
- (c) if a director signifies his agreement to the proposed written resolution, an alternate director appointed by him need not also signify his agreement; and
- (d) if an alternate director signifies his agreement to the proposed written resolution, his appointor need not also signify his agreement.

115 Proceedings of committees

115.1 All committees of the board shall, in the exercise of the powers delegated to them and in the transaction of business, conform with any mode of proceedings and regulations which the board may prescribe and subject thereto shall be governed by such of these articles as regulate the proceedings of the board as are capable of applying.

116 Minutes of proceedings

116.1 The board shall cause minutes to be made in books kept for the purpose of recording:

- (a) all appointments of officers and committees made by the board; and
- (b) the names of directors present at every meeting of the board, of a committee of the board, of the company or of the holders of any class of shares or debentures of the company, and all orders, resolutions and proceedings of such meetings.

116.2 Any such minutes, if purporting to be signed by the chairman of the meeting at which the proceedings were held or by the chairman of the next succeeding meeting or the secretary, shall be *prima facie* evidence of the matters stated in such minutes without any further proof.

117 Validity of proceedings

- 117.1 All acts done by a meeting of the board, or of a committee of the board, or by any person acting as a director, alternate director or member of a committee shall, notwithstanding that it is afterwards discovered that there was some defect in the appointment of any person or persons acting as aforesaid, or that they or any of them were or was disqualified from holding office or not entitled to vote, or had in any way vacated their or his office, be as valid as if every such person had been duly appointed, and was duly qualified and had continued to be a director, alternate director or member of a committee and entitled to vote.

DIRECTORS' INTERESTS

118 Director may have interests

- 118.1 A director may hold any other office or place of profit under the company (other than the office of auditor) in conjunction with his office of director on such terms as to tenure of office, remuneration and otherwise as the directors may determine. Any director may act by himself or his firm in a professional capacity (other than that of auditor) for the company and he or his firm shall be entitled to remuneration for such professional services.
- 118.2 Save as provided by the statutes, no director or intending director shall be disqualified by his office from contracting with the company either with regard to his tenure of any other office or place of profit, or as seller, buyer or otherwise; nor shall any such contract, or any contract or arrangement entered into by or on behalf of the company in which any director is in any way, whether directly or indirectly, interested, be liable to be avoided, nor shall any director so contracting or being so interested be liable to account to the company for any profit realised by any such contract or arrangement, by reason of such director holding that office or of the fiduciary relationship thereby established.
- 118.3
- (a) If a situation arises in which a director has, or can have, an interest that conflicts, or possibly may conflict, with the interests of the company (a “**relevant situation**”) the directors may, for the purposes of section 175 CA 2006, resolve to authorise:
- (i) if a relevant situation arises from the appointment or proposed appointment of a person as a director of the company, the appointment of the director and the relevant situation, subject to any limits or conditions which the directors may determine;
- (ii) if the relevant situation arises in circumstances other than as set out in article 118.3(a)(i), the relevant situation and the continuing

performance by the director of his duties, subject to any limits or conditions which the directors may determine,

and any such authorisation will be subject only to any limits or conditions which the directors expressly impose.

- (b) The interested director, and any other director with a similar interest, cannot vote, or be counted in the quorum, on a resolution to authorise his interest under article 118.3(a).
- (c) Any reference in article 118.3(a) to a conflict of interest includes a conflict of interest and duty and a conflict of duties.
- (d) Any limits or conditions determined by the directors under article 118.3(a) may be imposed at the time of authorisation or may be imposed or varied subsequently and may include (without limitation):
 - (i) whether the interested director(s) may vote (or be counted in the quorum at a meeting) in relation to any resolution relating to the relevant situation;
 - (ii) the exclusion of the interested director(s) from all information and discussion by the company of the relevant situation; and
 - (iii) the imposition of a specific duty of confidentiality for any confidential information of the company relating to the Relevant Situation.
- (e) An interested director must act in accordance with any limits or obligations imposed by the directors under article 118.3(a).
- (f) Subject to article 118.3(b), any authorisation under article 118.3(a) shall be dealt with in the same way as any other matter that may be decided by the directors under these articles.
- (g) Any authorisation of a relevant situation given by the directors under article 118.3(a) may provide that, where the interested director obtains (other than through his position as a director of the company) information that is confidential to a third party, he will not be obliged to disclose it to the company or to use it in relation to the company's affairs in circumstances where to do so would amount to a breach of that confidence.
- (h) Whilst there is a relevant situation, the general duties which the interested director owes to the company under sections 171 to 177 CA 2006 will not be infringed if he:

- (i) absents himself from meetings of the directors or from the discussion of any matter at a meeting relating to the relevant situation; and/or
- (ii) makes arrangements for papers to be received and read by a professional adviser on his behalf which may relate to the relevant situation; and/or
- (iii) behaves in any other way authorised by any guidance which may be issued by the directors from time to time.

118.4 (a) Any director may continue to be or become a director or other officer or member of or otherwise interested, as a member or otherwise, in any other company promoted by the company or any subsidiary undertaking of the company or in which the company or any subsidiary undertaking of the company may be interested or has any power of appointment and where such interest constitutes a relevant situation, such relevant situation shall be deemed to have been authorised pursuant to section 175 CA 2006, subject to any restrictions or conditions that may be imposed by the directors from time to time.

- (b) No such director shall be accountable for any remuneration or other benefits whatsoever received by him as a director or other officer or member of or from his interest in any such other company.
- (c) The directors may exercise the voting power conferred by the shares in any other company held or owned by the company, or exercisable by them as directors of such other company, in such manner in all respects as they think fit but subject to the like restrictions as are contained in article 118.6.

118.5 A director who is in any way, whether directly or indirectly, interested or deemed by the statutes to be interested in a contract, transaction or arrangement or a proposed contract, transaction or arrangement with the company shall declare the nature of his interest at a meeting of the directors in accordance with the statutes.

118.6 Save as provided in these articles and subject to the terms of any authorisation pursuant to article 118.3(a), a director shall not vote on or in respect of any contract or arrangement or any other proposal in which he has any interest which is to his knowledge a material interest otherwise than by virtue of his interests in shares or debentures or other securities or rights of or otherwise in or through the company. However a director shall be entitled to vote in respect of any contract or arrangement or any other proposal in which he has any interest which is not material. A director shall not be counted in the quorum at a meeting in relation to any resolution on which he is debarred from voting.

118.7 Subject to the provisions of the statutes a director (including an alternate director) shall (in the absence of some other material interest than is indicated below) be entitled to vote (and be counted in the quorum) in respect of a resolution relating to any of the following matters, namely:

- (a) the giving of any security, guarantee or indemnity in respect of:
 - (i) money lent or obligations incurred by him or any other person at the request of or for the benefit of the company or any of its subsidiary undertakings; or
 - (ii) a debt or obligation of the company or any of its subsidiary undertakings for which he himself has assumed responsibility in whole or in part under a guarantee or indemnity or by the giving of security; or
- (b) any proposal whereby the company or any of its subsidiary undertakings is offering securities under an offer in which he is or may be entitled to participate as a holder of securities or in the underwriting or sub-underwriting of which he is to participate; or
- (c) any proposal relating to any other company in which he does not to his knowledge hold an interest in shares (as that expression is defined for the purposes of Part VI of CA 2006) representing one per cent or more of either any class of the equity share capital of such company or the voting rights in such company; or
- (d) any arrangement for the benefit of employees of the company or any of its subsidiary undertakings which does not award him any privilege or benefit not generally awarded to the employees to whom such arrangement relates; and
- (e) any proposal concerning insurance which the company proposes to maintain or purchase for the benefit of persons including directors of the company.

118.8 Where proposals are under consideration concerning the appointment (including determining or varying the terms of appointment) of two or more directors to offices or employments with the company or any company in which the company is interested, such proposals may be divided and considered in relation to each director separately, and in such case each of the directors concerned shall, if not debarred from voting because of the limit on shareholding specified in article 118.7(c), be entitled to vote (and be counted in the quorum) in respect of each resolution except that concerning his own appointment.

- 118.9 If any question shall arise at any time as to the materiality of a director's interest or as to the entitlement of any director to vote and such question is not resolved by his voluntarily agreeing to abstain from voting, such question shall (subject to the statutes) be referred to the chairman of the meeting (or, where such question shall arise concerning such chairman, to such other director present at the meeting as the directors present, other than such chairman, shall by majority vote appoint) and his ruling in relation to any other director shall be final and conclusive except in a case where the nature or extent of the interests of the director concerned has not been fairly disclosed.
- 118.10 For the purposes of this article 118 the interest of any person who is connected with a director (within the meaning of section 252 CA 2006) shall be taken to be the interest of that director;
- 118.11 The directors may establish, maintain, participate in or contribute to, or procure the establishment and maintenance of, participation in or contribution to, any pension, annuity, superannuation, benevolent or life assurance fund, scheme or arrangement (whether contributory or otherwise) for the benefit of, and give or procure the giving of donations, gratuities, pensions, allowances, benefits and emoluments to, any persons who are or were at any time in the employment or service of the company, or any of its predecessors in business, or of any company which is a subsidiary undertaking of the company or is allied to or associated with the company, or with any such subsidiary undertaking, or who may be or have been directors or officers of the company, or of any such other company as previously stated, and the spouses, former spouses, families and dependants of any such persons, and also establish, subsidise and subscribe to any institutions, associations, societies, clubs, trusts or funds calculated to be for the benefit of or to advance the interests and well-being of the company or of any such other company as previously stated, or of any such persons as previously stated, and make payments for or towards the insurance of any such persons, and (subject to the provisions of the statutes) establish and contribute to any scheme for the acquisition of shares in the company or its holding company (whether or not an employees' share scheme) and (subject as previously stated) lend money to the company's employees to enable them to acquire such shares, and subscribe or guarantee money for charitable or benevolent objects, or for any exhibition or for any public, general or useful object, and do any of the matters previously stated either alone or in conjunction with others. Subject always, if the statutes shall so require, to particulars with respect to the proposed payment being disclosed to the members of the company and to the proposal being approved by ordinary resolution, any director shall be entitled to participate in and retain for his own benefit any such donation, gratuity, pension, allowance, benefit or emolument.

SEALS

119 Safe custody

119.1 The board shall provide for the safe custody of the seal, the securities seal and of any other seal of the company.

120 Application of seals

120.1 The seal shall be used only by the authority of a resolution of the board or of a committee of the board so authorised, which resolution may include a resolution for the giving of a general authority to any director, the secretary, an assistant secretary or other officer of the company to affix the seal from time to time. The board may determine whether any instrument to which the seal is affixed shall be signed and, if it is to be signed, who shall sign it and by what means. The board may also determine, either generally or in a particular case, that a signature may be dispensed with or affixed by mechanical, electronic or other means.

120.2 Every certificate or share warrant shall be issued either:

- (a) by affixing the securities seal to it, by mechanical, electronic or other means;
- (b) by printing a representation of the securities seal on it, by mechanical, electronic or other means, including laser printing; or
- (c) in such other manner as the board, having regard to the statutes and the regulations governing the Alternative Investment Market of the London Stock Exchange may authorise.

120.3 An instrument executed by a director and the secretary or by two directors with authority of a resolution of the board shall have the same effect as if executed under the seal.

121 Official seal for use abroad

121.1 Subject to the provisions of the statutes, the company may have an official seal for use in any place abroad.

THE SECRETARY

122 The secretary

122.1 Subject to the provisions of the statutes, the board shall appoint a secretary or joint secretaries and shall have power to appoint one or more persons to be an assistant or deputy secretary at such remuneration and on such terms and conditions as it thinks fit and any such person so appointed may be removed by the board. Such removal shall be without prejudice to any claim which any such

person may have for damages for breach of any contract of service between such person and the company.

- 122.2 Any provision of the statutes or of these articles requiring or authorising a thing to be done by or to a director and the secretary shall not be satisfied by its being done by or to the same person acting both as director and as the secretary.

AUTHENTICATION OF DOCUMENTS

123 Power to authenticate

- 123.1 Any director or the secretary or any person appointed for the purpose shall have power to authenticate any documents affecting the constitution of the company and any resolutions passed by the company or the directors or any committee of the directors, and any books, records, documents and accounts relating to the business of the company, and to certify copies thereof or extracts therefrom as true copies or extracts. Where any books, records, documents or accounts are elsewhere than at the office the local manager or other officer of the company having the custody thereof shall be deemed to be a person appointed by the directors as aforesaid. A document purporting to be a copy of a resolution, or an extract from the minutes of a meeting, of the company or of the directors or any committee of the directors which is certified as aforesaid shall be conclusive evidence in favour of all persons dealing with the company upon the faith thereof that such resolution has been duly passed or, as the case may be, that such minutes or extract is a true and accurate record of proceedings at a duly constituted meeting.

DIVIDENDS AND OTHER PAYMENTS

124 Declaration of dividends

- 124.1 Subject to the provisions of the statutes and of these articles, the company may by ordinary resolution declare dividends to be paid to members according to their respective rights and interests in the profits of the company. However, no dividend shall exceed the amount recommended by the board.

125 Interim dividends

- 125.1 Subject to the provisions of the statutes, the board may declare and pay such interim dividends (including any dividend payable at a fixed rate) as appears to the board to be justified by the profits of the company available for distribution. If at any time the share capital of the company is divided into different classes, the board may pay such interim dividends on shares which rank after shares conferring preferential rights with regard to dividend as well as on shares conferring preferential rights, unless at the time of payment any preferential dividend is in arrear. Provided that the board acts in good faith, it shall not incur any liability to the holders of shares conferring preferential rights for any loss that

they may suffer by the lawful payment of any interim dividend on any shares ranking after those with preferential rights.

126 Entitlement to dividends

126.1 Except as otherwise provided by the rights attached to shares, all dividends shall be declared and paid according to the amounts paid up (otherwise than in advance of calls) on the shares on which the dividend is paid. Subject as aforesaid, all dividends shall be apportioned and paid proportionately to the amounts paid up on the shares during any portion or portions of the period in respect of which the dividend is paid, but if any share is issued on terms providing that it shall rank for dividend as from a particular date, it shall rank for dividend accordingly.

127 Calls or debts may be deducted from dividends

127.1 The board may deduct from any dividend or other money payable to any person on or in respect of a share all such sums as may be due from him to the company on account of calls or otherwise in relation to the shares of the company.

128 Distribution in specie

128.1 The board may, with the authority of an ordinary resolution of the company, direct that payment of any dividend declared may be satisfied wholly or partly by the distribution of assets, and in particular of paid up shares or debentures of any other company, or in any one or more of such ways. Where any difficulty arises in regard to such distribution, the board may settle it as it thinks fit. In particular, the board may:

- (a) issue fractional certificates (or ignore fractions);
- (b) fix the value for distribution of such assets or any part thereof and determine that cash payments may be made to any members on the footing of the value so fixed, in order to adjust the rights of members; and
- (c) vest any such assets in trustees on trust for the persons entitled to the dividend.

129 Dividends not to bear interest

129.1 Unless otherwise provided by the rights attached to the share, no dividend or other moneys payable by the company or in respect of a share shall bear interest as against the company.

130 Method of payment

130.1 The company may pay any dividend, interest or other sum payable in respect of a share in cash or by direct debit, bank transfer, cheque, dividend warrant or money

order, by any method provided by the rules of a relevant system, or by any other method (including by electronic media) as the board may consider appropriate and may send the same by post or other delivery service (or by such other means offered by the company as the member or persons entitled to it may agree in writing) to the registered address (or in the case of a depository, subject to the approval of the board, such persons and addresses as the depository may require) of the member or person entitled to it (or, if two or more persons are holders of the share or are jointly entitled to it by reason of the death or bankruptcy of the member or otherwise by operation of law, to the registered address of such of those persons as is first named in the register) or to such person and such address as such member or person or persons may direct in writing.

130.2 Every cheque, warrant, order or other form of payment is sent at the risk of the person entitled to the money represented by it, shall (where relevant) be crossed in accordance with the Cheques Act 1992 and shall be made payable to the person or persons entitled, or to such other person as the person or persons entitled may direct in writing. Payment of the cheque, warrant, order or other form of payment shall be a good discharge to the company. If any such cheque, warrant, order or other form of payment has or shall be alleged to have been lost, stolen or destroyed, the board may, at the request of the person entitled thereto, issue a replacement cheque or warrant or order or make payment in some other form, subject to compliance with such conditions as to evidence and indemnity and the payment of out of pocket expenses of the company in connection with the request as the board may think fit.

130.3 Any joint holder or other person jointly entitled to a share may give an effective receipt for any dividend or other moneys payable in respect of such share.

130.4 The board may, at its discretion, make provisions to enable a depository and/or any member as the board shall from time to time determine to receive duly declared dividends in a currency or currencies other than sterling. For the purposes of the calculation of the amount receivable in respect of any dividend, the rate of exchange to be used to determine the foreign currency equivalent of any sum payable as a dividend shall be such rate or rates and the payment thereof shall be on such terms and conditions as the board may in its absolute discretion determine.

131 Uncashed dividends

131.1 If cheques, warrants or orders for dividends or other sums payable in respect of a share sent by the company to the person entitled thereto are returned to the company or left uncashed on two consecutive occasions or, following one occasion, reasonable enquiries have failed to establish any new address to be used for the purpose, the company shall not be obliged to send any dividends or other moneys payable in respect of that share due to that person until he notifies the company of an address to be used for the purpose.

132 Unclaimed dividends

132.1 All dividends, interest or other sum payable and unclaimed for twelve months after having become payable may be invested or otherwise made use of by the board for the benefit of the company until claimed and the company shall not be constituted a trustee in respect thereof. All dividends unclaimed for a period of twelve years after having been declared or become due for payment shall (if the board so resolves) be forfeited and shall cease to remain owing by the company.

133 Payment of share dividends

133.1 The board may, with the prior authority of an ordinary resolution of the company and subject to such terms and conditions as the board may determine, offer to any holders of ordinary shares the right to elect to receive ordinary shares, credited as fully paid, instead of the whole (or some part, to be determined by the board) of any dividend specified by the ordinary resolution. The following provisions shall apply:

- (a) the said resolution may specify a particular dividend, or may specify all or any dividends declared within a specified period or periods;
- (b) the entitlement of each holder of ordinary shares to new ordinary shares shall be such that the relevant value of the entitlement shall be as nearly as possible equal to (but not greater than) the cash amount (disregarding any tax credit) of the dividend that such holder would have received by way of dividend. For this purpose "relevant value" shall be calculated by reference to the average of the middle market quotations for the ordinary shares on the Alternative Investment Market of the London Stock Exchange for the day on which the ordinary shares are first quoted "ex" the relevant dividend and the four subsequent dealing days, or in such other manner as may be determined by or in accordance with the ordinary resolution. A certificate or report by the auditors as to the amount of the relevant value in respect of any dividend shall be conclusive evidence of that amount;
- (c) no fractions of a share shall be allotted. The board may make such provisions as it thinks fit for any fractional entitlements, including provisions whereby, in whole or in part, the benefit thereof accrues to the company and/or under which fractional entitlements are accrued and/or retained and in each case accumulated on behalf of any member and such accruals or retentions are applied to the allotment by way of bonus to or cash subscription on behalf of such member of fully paid ordinary shares and/or provisions whereby cash payments may be made to members in respect of their fractional entitlements;

- (d) the board shall, after determining the basis of allotment, notify the holders of ordinary shares in writing of the right of election offered to them, and specify the procedure to be followed and place at which, and the latest time by which, elections must be lodged in order to be effective;
- (e) the board may exclude from any offer any holders of ordinary shares or any ordinary shares held by a depositary or any ordinary shares on which dividends are payable in foreign currency where the board considers that the making of the offer to them or in respect of such shares would or might involve the contravention of the laws of any territory or that for any other reason the offer should not be made to them or in respect of such shares;
- (f) the board may establish or vary from time to time a procedure for election mandates in respect of future rights of election and may determine that every duly effected election in respect of any ordinary shares shall be binding on every successor in title to the holder thereof;
- (g) the dividend (or that part of the dividend in respect of which a right of election has been offered) shall not be payable on ordinary shares in respect of which an election has been duly made (the “**elected ordinary shares**”) and instead additional ordinary shares shall be allotted to the holders of the elected ordinary shares on the basis of allotment determined as aforesaid. For such purpose the board may capitalise, out of any amount for the time being standing to the credit of any reserve or fund (including any share premium account, capital redemption reserve or other undistributable reserve) or of any of the profits which could otherwise have been applied in paying dividends in cash as the board may determine, a sum equal to the aggregate nominal amount of the additional ordinary shares to be allotted on that basis and apply it in paying up in full the appropriate number of unissued ordinary shares for allotment and distribution to the holders of the elected shares on that basis. A board resolution capitalising any part of such reserve or fund or profits shall have the same effect as if such capitalisation had been declared by ordinary resolution of the company in accordance with article 135 and in relation to any such capitalisation the board may exercise all powers conferred on it by article 135 without need of such ordinary resolution;
- (h) the additional ordinary shares so allotted shall rank *pari passu* in all respects with each other and with the fully paid ordinary shares in issue on the record date for the dividend in respect of which the right of election has been offered, except that they will not rank for any dividend or other distribution or other entitlement which has been declared, paid or made by reference to such record date; and
- (i) the board may terminate, suspend or amend any offer of the right to elect to receive ordinary shares in lieu of any cash dividend at any time and

generally may implement any share dividend scheme on such terms and conditions as the board may from time to time determine and take such other action as the board may deem necessary or desirable from time to time in respect of any such scheme.

134 Reserves

134.1 The board may, before recommending any dividend (whether preferential or otherwise), carry to reserve out of the profits of the company such sums as it thinks fit. All sums standing to reserve may be applied from time to time, at the discretion of the board, for any purpose to which the profits of the company may properly be applied, and pending such application may, at the like discretion, either be employed in the business of the company or be invested in such investments as the board thinks fit. The board may divide the reserve into such special funds as it thinks fit and may consolidate into one fund any special funds or any parts of any special funds into which the reserve may have been divided as it thinks fit. Any sum which the board may carry to reserve out of the unrealised profits of the company shall not be mixed with any reserve to which profits available for distribution have been carried. The board may also, without placing the same to reserve, carry forward any profits which it may think prudent not to distribute.

135 Capitalisation of reserves

135.1 The board may, with the authority of an ordinary resolution of the company:

- (a) subject as provided in this article, resolve to capitalise any undivided profits of the company not required for paying any preferential dividend (whether or not they are available for distribution) or any sum standing to the credit of any reserve or fund of the company which is available for distribution or standing to the credit of share premium account or capital redemption reserve or other undistributable reserve;
- (b) appropriate the sum resolved to be capitalised to the holders of ordinary shares in proportion to the nominal amounts of the shares (whether or not fully paid) held by them respectively which would entitle them to participate in a distribution of that sum if the shares were fully paid and the sum were then distributable and were distributed by way of dividend and apply such sum on their behalf either in or towards paying up the amounts, if any, for the time being unpaid on any shares held by them respectively, or in paying up in full unissued shares or debentures of the company of a nominal amount equal to that sum, and allot the shares or debentures credited as fully paid to those holders of ordinary shares or as they may direct, in those proportions, or partly in one way and partly in the other, provided that:

- (i) the share premium account, the capital redemption reserve, any other undistributable reserve and any profits which are not available for distribution may, for the purposes of this article, only be applied in paying up unissued shares to be allotted to holders of ordinary shares credited as fully paid; and
 - (ii) in a case where any sum is applied in paying amounts for the time being unpaid on any shares of the company or in paying up in full debentures of the company, the amount of the net assets of the company at that time is not less than the aggregate of the called up share capital of the company and its undistributable reserves as shown in the latest audited accounts of the company or such other accounts as may be relevant and would not be reduced below that aggregate by the payment thereof;
- (c) resolve that any shares so allotted to any member in respect of a holding by him of any partly paid shares shall, so long as such shares remain partly paid, rank for dividends only to the extent that such partly paid shares rank for dividends;
- (d) make such provision by the issue of fractional certificates (or by ignoring fractions or by accruing the benefit thereof to the company rather than to the holders of ordinary shares concerned) or by payment in cash or otherwise as it thinks fit in the case of shares or debentures becoming distributable in fractions;
- (e) authorise any person to enter on behalf of all the holders of ordinary shares concerned into an agreement with the company providing for either:
 - (i) the allotment to them respectively, credited as fully paid up, of any shares or debentures to which they may be entitled on such capitalisation; or
 - (ii) the payment up by the company on behalf of such holders by the application thereto of their respective proportions of the reserves or profits resolved to be capitalised, of the amounts or any part of the amounts remaining unpaid on their existing shares,any agreement made under such authority being effective and binding on all such holders); and
- (f) generally do all acts and things required to give effect to such resolution.

136 Record dates

136.1 Notwithstanding any other provision of these articles but without prejudice to the rights attached to any shares and subject always to the statutes, the company or the board may by resolution specify any date (the "**record date**") as the date at the close of business (or such other time as the board may determine) on which persons registered as the holders of shares or other securities shall be entitled to receipt of any dividend, distribution, interest, allotment, issue, notice, information, document or circular and such record date may be on or at any time before the date on which the same is paid, made, given or served or (in the case of any dividend, distribution, interest, allotment or issue) at any time after the same is recommended, resolved, declared or announced, but without prejudice to the rights *inter se* in respect of the same of transferors and transferees of any such shares or other securities. No change in the register of such holders after the record date shall invalidate the same.

ACCOUNTS

137 Inspection of records

137.1 No member (other than a director) shall have any right to inspect any accounting record or other document of the company unless he is authorised to do so by statute, by order of the court, by the board or by ordinary resolution of the company.

138 Accounts to be sent to members

138.1 Except as provided in article 139, a copy of the directors' and auditors' reports accompanied by copies of the annual accounts shall, not less than twenty one clear days before the annual general meeting before which they are to be laid, be delivered or sent or made available to every member and holder of debentures of the company and to the auditors and to every other person who is entitled to receive notice of general meetings. If and to the extent permitted by the statutes, all or any of such documents may be delivered to a member by means of electronic mail. However, this article shall not require a copy of those documents to be sent or made available to any person who, under the provisions of these articles, is not entitled to receive notices from the company or of whose address the company is unaware or to any holder of debentures of whose address the company is unaware or to more than one of the joint holders of any shares or debentures. If all or any of the shares in or debentures of the company are listed or dealt in on any stock exchange, there shall at the same time be forwarded to the secretary of that stock exchange such number of copies of each of those documents as the regulations of that stock exchange may require.

139 Summary financial statements

139.1 The company may, in accordance with section 251 CA 1985 and any regulations made under it, send a summary financial statement to any member instead of or in addition to the documents referred to in article 138. Where it does so, the statement shall be delivered or sent by post to the member not less than twenty one clear days before the annual general meeting before which those documents are to be laid. If and to the extent permitted by the statutes, such summary financial statement may be delivered by means of electronic mail.

NOTICES

140 Notices to be in writing

140.1 Any notice to be given to or by any person pursuant to these articles shall be in writing, except that a notice convening a board or board committee meeting need not be in writing.

141 Service of notice on members

141.1 Subject to article 140.1 and unless otherwise provided by these articles, the company shall send or supply a document or information that is required or authorised to be sent or supplied to a member or any other person by the company by a provision of the statutes or pursuant to these articles or by any other rules or regulations to which the company may be subject, in such form and by such means as the board may in its absolute discretion determine, provided that the provisions of CA 2006 which apply to a document or information required or authorised to be sent or supplied by the statutes shall, in the same way, also apply to any document or information required or authorised to be sent by these articles or by any other rules or regulations to which the company may be subject.

141.2 Subject to article 140.1 and unless otherwise provided by these articles, a member or a person entitled by transmission to a share shall send a document or information to the company pursuant to these articles in such form and by such means as it may in its absolute discretion decide, provided that the determined form and means are permitted by the statutes for the purpose of sending or supplying a document or information of that type to a company pursuant to a provision of the statutes and, unless the board otherwise permits, any applicable condition or limitation specified by the statutes, including without limitation as to the address to which the document or information may be sent, is satisfied and, unless otherwise provided by these articles or required by the board, such document or information shall be authenticated in the manner specified by the statutes for authentication of a document or information sent in the relevant form.

- 141.3 In the case of joint holders of a share, all notices or documents shall be given to the joint holder whose name stands first in the register in respect of the joint holding. Notice so given shall be sufficient notice to all the joint holders.
- 141.4 Where a member (or, in the case of joint holders, the person first named in the register) has a registered address which is not within an EEA State but has notified the company of an address within an EEA State at which notices, documents or other information may be sent to him, he shall be entitled to have notices, documents or other information sent to him at that address (provided, in the case of a notice, document or information sent by electronic means, including without limitation any notification required by the statutes that the document or information is available on a website, the company so agrees, which agreement the company shall be entitled in its absolute discretion to withhold, including without limitation in circumstances in which the company considers that the sending of the document or information to such address using electronic means would or might infringe the laws of any jurisdiction); but otherwise no such member shall be entitled to receive any notice, document or information from the company and, without prejudice to the generality of the foregoing, any notice of an annual general meeting or general meeting which is in fact sent or purports to be sent to such member shall be ignored for the purpose of determining the validity of the proceedings at such meeting.
- 141.5 If on three consecutive occasions notices or other documents have been sent through the post to any member at his registered address or his address for the service of notices but have been returned undelivered, such member shall not thereafter be entitled to receive notices or other documents from the company until he shall have communicated with the company and supplied in writing a new registered address or address within an EEA State for the service of notices.
- 141.6 If on two consecutive occasions notices or other documents have attempted to be sent in electronic form to any member at an address for the time being notified by that member to the company for such purpose and the company is aware that there has been a failure of delivery of such notice or document, then the company shall thereafter send notices or documents to that member through the post to his registered address or his address for service of notices by post, in which case the provisions of article 141.5 shall apply.
- 141.7 The board may from time to time issue, endorse or adopt terms and conditions relating to the use of electronic means for the sending of notices, documents or other information, including proxy appointments, by the company to members or persons entitled by transmission and by members or persons entitled by transmission to the company.
- 141.8 Without prejudice to articles 141.1 and 141.2 but subject to CA 2006, notices, documents and other information shall be validly sent or supplied by the company if they are made available on a website. Notices, documents or other information

sent to the company in electronic form shall not be treated as received by the company if they are rejected by computer virus protection arrangements.

142 Notice in case of Death, Bankruptcy or Mental Disorder

142.1 The company may give notice to the person entitled to a share in consequence of the death or bankruptcy of a member or otherwise by operation of law, by sending or delivering it in any manner authorised by these articles for the giving of notice to a member, addressed to that person by name, or by the title of representative of the deceased or trustee of the bankrupt or representative by operation of law or by any like description, at the address (if any) within the United Kingdom supplied for the purpose by the person claiming to be so entitled. Until such an address has been so supplied, a notice may be given in any manner in which it might have been given if the death or bankruptcy or operation of law had not occurred.

143 Evidence of Service

143.1 Any member present, in person or by proxy, at any meeting of the company or of the holders of any class of shares of the company shall be deemed to have received due notice of such meeting, and, where requisite, of the purposes for which such meeting was called.

143.2 Proof that a notice, document or other information was properly addressed, prepaid and posted shall be conclusive evidence that the notice, document or information was sent.

143.3 A notice, document or information sent by the company to a member by post shall be deemed to have been received:

(a) if sent by first class post or special delivery post from an address in the United Kingdom to another address within the United Kingdom, or by a postal service similar to first class post or special delivery post from an address in another country to another address in that same other country, on the day following that on which the notice, document or information was posted;

(b) if sent by airmail from an address in the United Kingdom to an address outside the United Kingdom, or from an address in another country to an address outside that country (including without limitation an address within the United Kingdom), on the third day following that on which the notice, document or information was posted; and

(c) in any other case, on the second day following that on which the notice, document or information was posted.

143.4 Proof that a notice, document or other information sent or supplied by electronic means was properly addressed and despatched shall be conclusive evidence that

the notice, document or other information was sent or supplied, unless the company is aware that there has been a failure of delivery of such notice, document or information following two attempts, in which case such notice, document or information shall be sent to the member at his registered address or address for service, provided that the date of deemed service or delivery shall be 48 hours after the time that such notice, document or information was originally sent by electronic means in accordance with article 143.5.

143.5 A notice, document or information sent or supplied by the company to a member in electronic form shall be deemed to have been received by that member at the expiration of 48 hours after the time that such notice, document or information was sent or supplied to that member. Such notice, document or information shall be deemed received by the member on that day notwithstanding that the company becomes aware that the member has failed to receive the relevant document or information for any reason and notwithstanding that the company subsequently sends a hard copy of such notice, document or information by post to the member.

143.6 A notice, document or information sent or supplied by the company to a member by means of a website shall be deemed to have been received by that member:

(a) when the notice, document or information was first made available on the website; or

(b) if later, when the member is deemed by article 143.5 or article 143.3 to have received notice of the fact that the notice, document or information was available on the website; and such a notice, document or information shall be deemed received by the member on that day notwithstanding that the company becomes aware that the member has failed to receive the relevant notice, document or information for any reason and notwithstanding that the company subsequently sends a hard copy of such notice, document or information by post to the member.

144 Notice Binding on Transferees

144.1 Every person who, by operation of law, transfers or by any other means becomes entitled to a share shall be bound by any notice in respect of that share (other than a section 793 notice) which, before his name is entered in the register, has been duly given to a person from whom he derives his title.

145 Notice by Advertisement

145.1 Any notice to be given by the company to the members or any of them, and not otherwise provided for by these articles, shall be sufficiently given if given by advertisement in at least one national newspaper published in the United Kingdom and, where the company keeps an overseas branch register, in at least one daily newspaper published in the territory in which such register is maintained. Any

notice given by advertisement shall be deemed to have been served at noon on the day on which the advertisement first appears.

146 Suspension of Postal Services

- 146.1 If at any time by reason of the suspension, interruption or curtailment of postal services or threat thereof within the United Kingdom the company is or would be unable effectively to convene an annual general meeting or a general meeting by notices sent through the post, such a meeting may be convened by a notice advertised in at least one national newspaper published in the United Kingdom and, where the company keeps an overseas branch register, in at least one daily newspaper published in the territory in which such register is maintained. Such notice shall be deemed to have been duly served on all members entitled thereto at noon on the day on which the first of such advertisements appears. In any such case the company shall send confirmatory copies of the notice by post to any member who has not previously been sent the notice by another method permitted by these articles if, at least seven days prior to the meeting, the posting of notices to addresses throughout the United Kingdom again becomes practicable.

WINDING UP

147 Division of assets

- 147.1 If the company is wound up the liquidator may, with the sanction of a special resolution of the company and any other sanction required by law, divide among the members in specie the whole or any part of the assets of the company and may, for that purpose, value any assets and determine how the division shall be carried out as between the members or different classes of members. Any such division may be otherwise than in accordance with the existing rights of the members, but if any division is resolved otherwise than in accordance with such rights, the members shall have the same right of dissent and consequential rights as if such resolution were a special resolution passed pursuant to section 110 of the Insolvency Act 1986. The liquidator may, with the like sanction, vest the whole or any part of the assets in trustees on such trusts for the benefit of the members as he with the like sanction shall determine, but no member shall be compelled to accept any assets on which there is a liability.

148 Transfer or sale under section 110 Insolvency Act 1986

- 148.1 A special resolution sanctioning a transfer or sale to another company duly passed pursuant to section 110 of the Insolvency Act 1986 may in the like manner authorise the distribution of any shares or other consideration receivable by the liquidator among the members otherwise than in accordance with their existing rights, and any such determination shall be binding on all the members, subject to the right of dissent and consequential rights conferred by the said section.

INDEMNITY

149 Right to indemnity

149.1 Without prejudice to any indemnity to which he may otherwise be entitled (and subject to article 149.4), every person who is or was, at any time on or after 1 January 2006, a director, secretary or other relevant officer of the company shall be indemnified and kept indemnified out of the company's assets against all liability incurred by him as such or as a director, secretary or relevant officer of an associated company, including where such company is trustee of an occupational pension scheme:

- (a) in defending any proceedings, whether civil or criminal, in respect of alleged negligence, default, breach of duty, breach of trust or otherwise in relation to the company or an associated company or its or their affairs, in which judgement is given in his favour or in which he is acquitted or in defending or settling any such proceedings which are otherwise disposed of on terms previously agreed with the board or on terms otherwise approved by the board without a finding or admission of negligence, default, breach of duty or breach of trust on his part; or
- (b) in connection with any application under the statutes in which relief is granted to him by the court or any investigations or actions of a regulatory authority as referred to in section 206 CA 2006,

provided that this article shall not grant, or entitle any such person to, indemnification to the extent that it would cause this article, or any part of it, to be void under the statutes.

149.2 Without prejudice to any indemnity to which he may otherwise be entitled (including, for the avoidance of doubt, any indemnity under or pursuant to these articles) and to the extent permitted by the statutes, the board shall have power in the name and on behalf of the company to:

- (a) grant on such terms as it sees fit any person who is or was, at any time on or after 1 January 2006, a director, secretary or other relevant officer of the company an indemnity or indemnities out of the assets of the company in respect of any liability incurred by him as such or as a director, secretary or relevant officer of an associated company, including where such company is a trustee of an occupational pension scheme, and to amend, vary or extend the terms of any such indemnity so granted, again on such terms as the board sees fit; and/or
- (b) enter into and amend, vary or extend such arrangements as it sees fit:
 - (i) to provide any person who is or was, at any time on or after 1 January 2006, a director, secretary or other relevant officer of the

company with funds to meet expenditure incurred or to be incurred by him in defending any criminal or civil proceedings brought against him as such or as a director, secretary or relevant officer of an associated company or in connection with any application for relief under the statutes or any investigation or actions of a regulatory authority as referred to in section 206 CA 2006; or

- (ii) to enable any such person to avoid incurring any such expenditure.

149.3 For the purposes of this article 149:

- (a) a "**relevant officer**" is any officer of the company or an associated company (other than in either case any person (whether or not an officer of the company or an associated company) engaged by the company or an associated company as auditor);
- (b) "**associated company**" has the meaning given to that term in section 256 CA 2006; and
- (c) a director shall be entitled to vote and to be counted in the quorum at any meeting of the board or a committee of the board at which any indemnity, arrangement or proposal falling within any of the provisions of articles 149.1 or 149.2 is to be considered and, for the purposes of article 118, any interest which any director may have in such indemnity, arrangement or proposal shall not be a material interest unless the terms of such indemnity arrangement or proposal confer upon such director a privilege or benefit not generally available to, or awarded to, any other director. The decision of the chairman of the meeting as to whether the indemnity, arrangement or proposal to be considered at the meeting falls within the provisions of articles 149.1 or 149.2 or as to the materiality of any director's interest therein for the purposes of this article and article 118 shall be final and conclusive.

149.4 Subject to the provisions of the statutes but without prejudice to any indemnity to which he may otherwise be entitled, every person who ceased to be a director or other officer or auditor of the company on or before 31 December 2005 shall be indemnified out of the assets of the company against any liability incurred by him in defending any proceedings, whether civil or criminal, in which judgment is given in his favour or in which he is acquitted or in connection with any application in which relief is granted to him by the court from liability for negligence, default, breach of duty or breach of trust in relation to the affairs of the company.