



Cyprotex PLC,
15 Beech Lane,
Macclesfield,
Cheshire SK10 2DR
Tel: +44 (0)1625 505 100
Fax: +44 (0)1625 505 199

19th June 2008

Dear Shareholder

Annual General Meeting: 14 July 2008

I am pleased to enclose the annual report and accounts for your Company to 31 December 2007 together with the notice of the annual general meeting (AGM) and a proxy card.

The resolutions to be proposed at the AGM relate to normal business except for two of the resolutions. As you may have seen from an announcement issued to the Stock Exchange on 3 June 2008, we have agreed terms with Robert Morrisson Atwater for the termination of his employment with the Group. At the same time we have agreed terms with Dr Anthony (Tony) Baxter for his appointment as our new CEO.

In accordance with company law, Tony Baxter's appointment needs to be confirmed at the AGM given that his appointment was made by the board.

Also, as required by company law, the approval of the shareholders to the ex gratia element of the compensation payable to Robert for his loss of office will be sought at the AGM. A memorandum summarising the terms of the settlement with Robert will be available for inspection at the Company's registered office 15 days before the date of the AGM and also at the meeting itself. If in the meantime you have any questions relating to Robert's settlement please let me know.

Pursuant to s.95 of the Companies Act 1985, the directors recommend resolution No.8 relating to the issue of up to £13,865 of shares free of pre-emption rights at such price and on such terms as they think fit. We believe that the flexibility given to your board in this respect is in the best interests of shareholders to help to try to drive the business forward.

I look forward to seeing you at the National Liberal Club at 1 Whitehall Place for the AGM on 14 July 2008 at 10.00am.

Yours faithfully

Nikolas Sofronis
Chairman, Cyprotex PLC