

THIS DOCUMENT IS IMPORTANT AND REQUIRES YOUR IMMEDIATE ATTENTION. Your attention is drawn to the Circular (as defined below). If you are in any doubt as to the action you should take, you should immediately consult your stockbroker, solicitor, accountant or other independent financial adviser authorised under the Financial Services and Markets Act 2000

CYPROTEX PLC

(the Company)

(Incorporated in England and Wales under the Companies Act 1985 with registered number 04311107)

APPLICATION FORM

Offer of new ordinary shares of 0.1p each ("New Ordinary Shares")

Terms defined in the Chairman's letter of 20 June 2008 and documents from the Company to the shareholders enclosed with this application form (the **Circular**) shall have the same meaning in this Application Form.

Box A	Your name and address <i>(please print)</i> <i>Name</i> <i>Address</i>
Box B	The number of Shares you hold as at the close of business on 11 July 2008 ("X") <i>Number</i>
Box C	The maximum number of New Ordinary Shares you are entitled to subscribe for is (using the following formula) ("Y"). Where: $Y = \left(\frac{178,647,988}{138,647,988} \times X \right) - X$ Or more simply $Y = 0.2885004 \times X$ <i>(your calculation of "Y" should be rounded up to the nearest whole number)</i> <i>Number</i>
Box D	Do you wish to subscribe for New Ordinary Shares? If so, for how many? ("Z")	Yes / No <i>(please delete as appropriate)</i> <i>Number (please round up to the nearest whole number)</i>

Box E	The amount payable for your entitlement Z x £0.025	£..... <i>Amount</i>
--------------	--	-------------------------

If you have any questions on the application procedures set out in this document, you should contact Russell Gibbs or Mark C. Warburton at the Company on 01625 505 100.

To: The Directors

I/We acknowledge receipt of the Circular and I/we confirm that, on the terms and subject to the conditions set out therein, I/we irrevocably subscribe for and accept the allocation to me/us of the number of shares stated in Box D subject to the terms set out in the Circular (**Placing Participation**).

I/We undertake, not later than midday on Tuesday, 26 August 2008 to make payment arranged either (i) by bank telegraphic transfer to Bank of Scotland, St. James’s Gate, 14/16 Cockspur Street, London SW1Y 5BL, account name: Cyprotex PLC, account number 06090664, sort code: 12-11-03, IBAN GB03 BOFS 1211 0306 0906 64, SWIFT (or BIC) BOFSGB21238 quoting “Shareholder Offer” or (ii) by cheque (payable to Cyprotex PLC), for the full amount due in respect of the number of New Ordinary Shares for which I/we have agreed to subscribe for at 2.5p per New Ordinary Share. If the conditions to which the Placing Participation is subject are not satisfied, I/we understand that my/our subscription monies will be returned to me/us (without interest) as soon as practicable thereafter.

I/we acknowledge that all documents and remittances relating to the Placing Participation are sent at my/our risk.

I/we hereby request you to procure the issue and delivery to me/us of a definitive certificate in respect of the Placing Participation in accordance with the instructions given below, at my/our risk.

I/We confirm that I/we have not relied on any information or representation relating to the Placing or the Company other than information which is contained in the Circular. By returning this Application Form I/we give all the warranties, undertakings, confirmations and acknowledgements specified in the Circular (and in particular the warranties, undertakings and confirmations set out in the paragraph headed “*Confirmations, warranties and undertakings*”).

I/we confirm and warrant that I am/we are not liable and I am/we are not accepting as nominee(s) or agent(s) for a person who is or may be liable to pay stamp duty at any rate higher than £5.00 for every £1,000 or part of £1,000 or stamp duty reserve tax at a rate higher than 0.5% of the 2.5p per New Ordinary Share which I/we will subscribe.

I/we also confirm that I/we have obtained all necessary consents and authorities to enable me/us to accept the Shareholder Offer and to perform my/our obligations specified in the Circular. I/We confirm that I/we will provide you with such information and documents as you

or the Company may require in order to comply with your obligations under the Money Laundering Regulations 2007 (as amended).

Completed Application Forms should be signed and dated and returned to the Company Secretary at the Company's office at 15 Beech Lane, Macclesfield, Cheshire SK10 2DR, together with a cheque (made payable to Cyprotex PLC) or telegraphic transfer to the bank account stated above, for payment in full for the appropriate amount (rounded up to the nearest whole pence), to be received by the Company IN CLEARED FUNDS by no later than midday on Tuesday, 26 August 2008.

Signed: *[Eligible Shareholder(s) should sign here]*

Date:2008

A corporation should execute this form acting in accordance with the laws of the territory in which the corporation is incorporated.