

# Cyprotex PLC

("Cyprotex" or the "Company")

Holding(s) in Company

TR-1: NOTIFICATION OF MAJOR INTEREST IN SHARES <sup>i</sup>	
<b>1. Identity of the issuer or the underlying issuer of existing shares to which voting rights are attached:</b> <sup>ii</sup>	Cyprotex PLC
<b>2 Reason for the notification</b> (please tick the appropriate box or boxes):	
An acquisition or disposal of voting rights	<input type="checkbox"/>
An acquisition or disposal of qualifying financial instruments which may result in the acquisition of shares already issued to which voting rights are attached	<input type="checkbox"/>
An acquisition or disposal of instruments with similar economic effect to qualifying financial instruments	<input type="checkbox"/>
An event changing the breakdown of voting rights	<input type="checkbox"/>
Other (please specify):	Transfer in to Nominees <input checked="" type="checkbox"/>
<b>3. Full name of person(s) subject to the notification obligation:</b> <sup>iii</sup>	Rensburg Sheppards Investment Management Limited
<b>4. Full name of shareholder(s)</b> (if different from 3.): <sup>iv</sup>	N/A
<b>5. Date of the transaction and date on which the threshold is crossed or reached:</b> <sup>v</sup>	17 <sup>th</sup> June 2010
6. Date on which issuer notified:	18 <sup>th</sup> June 2010
<b>7. Threshold(s) that is/are crossed or reached:</b> <sup>vi, vii</sup>	5%

8. Notified details:							
A: Voting rights attached to shares <sup>viii, ix</sup>							
Class/type of shares  if possible using the ISIN CODE	Situation previous to the triggering transaction		Resulting situation after the triggering transaction				
	Number of Shares	Number of Voting Rights	Number of shares	Number of voting rights		% of voting rights <sup>x</sup>	
			Direct	Direct <sup>xi</sup>	Indirect <sup>xii</sup>	Direct	Indirect
Ordinary Shares GB0031214470					10,401,600		5.81%
<b>B: Qualifying Financial Instruments</b>							
Resulting situation after the triggering transaction							

Type of financial instrument	Expiration date <sup>xiii</sup>	Exercise/Conversion Period <sup>xiv</sup>	Number of voting rights that may be acquired if the instrument is exercised/converted.	% of voting rights	
<b>C: Financial Instruments with similar economic effect to Qualifying Financial Instruments</b> <sup>xv, xvi</sup>					
Resulting situation after the triggering transaction					
Type of financial instrument	Exercise price	Expiration date <sup>xvii</sup>	Exercise/Conversion period <sup>xviii</sup>	Number of voting rights instrument refers to	% of voting rights <sup>xix, xx</sup>
					Nominal
					Delta
Total (A+B+C)					
Number of voting rights				Percentage of voting rights	
10,401,600				5.81%	

<b>9. Chain of controlled undertakings through which the voting rights and/or the financial instruments are effectively held, if applicable:</b> <sup>xxi</sup>
Client holdings registered in the name of Nominee companies 100% owned by Rensburg Sheppards Investment Management Limited.

Proxy Voting:	
10. Name of the proxy holder:	N/A
11. Number of voting rights proxy holder will cease to hold:	N/A
12. Date on which proxy holder will cease to hold voting rights:	N/A