

**THE COMPANIES ACTS 1985 AND 1989**

**PUBLIC COMPANY LIMITED BY SHARES**

**MEMORANDUM**

**OF**

**ASSOCIATION**

**OF**

**CYPROTEX PLC**



**ADDLESHAW BOOTH & CO**

Company No: 4311107

**THE COMPANIES ACTS 1985 AND 1989**

---

**PUBLIC COMPANY LIMITED BY SHARES**

---

**MEMORANDUM OF ASSOCIATION**

**OF**

**CYPROTEX PLC**

- 1 The Company's name is "Cyprotex PLC".<sup>1</sup>
- 2 The Company's registered office is to be situated in England and Wales.
- 3 The Company is a Public Limited Company.
- 4 The Company's objects are:
  - (a) To carry on the business of a holding company in all of its branches and to co-ordinate the policy and administration and deal with or assist in dealing with the administration and accounting functions of any companies of which the Company is a member or which are in any manner controlled by, or connected with, the Company and to establish or promote or concur in establishing or promoting any other company whose objects shall include the acquisition and taking over of all or any of the assets and liabilities of the Company or the promotion of which shall be in any manner calculated to advance directly or indirectly the objects or interests of the Company, and to acquire by purchase, lease, concession, grant, licence or otherwise and hold or dispose of such businesses, options, rights, privileges, lands, buildings, leases, underleases, debentures, bonds, obligations, reversionary interests, annuities, policies of assurance, shares, stock or securities and any other property as the Company shall deem fit, and guarantee with or without security the payment of dividends, interest or capital of any shares, stock or securities issued by, or any other obligations of, any such company.
  - (b) To carry on all or any of the businesses of general merchants and traders, cash and credit traders, importers, exporters, designers, manufacturers, manufacturers' agents and representatives, buyers, sellers, distributors, factors, wholesalers, retailers, letters on hire and shippers of and dealers in produce, products, goods,

---

<sup>1</sup> The Company's name on incorporation was Inhoco 2437 Limited. It was changed to Cyprotex Services Limited on 4 December 2001 and to Cyprotex Limited on 4 January 2002. On 4 January 2002 the Company was re-registered as a plc.

