

## Cerep welcomes President Clinton's proposals for the human genome

**Paris, France, 16 March 2000** – CEREP has welcomed the recent joint proposals by President Clinton and UK Prime Minister Tony Blair to keep information on the human genome in the public domain.

Unrestricted use of this information by the pharmaceuticals industry and biotechnology companies would be favourable for CEREP which, for over ten years, has specialised in streamlining pre-clinical trials in areas of pharmacological interest for drug discovery.

More generally, free access to the findings of genomic research will further enhance the development of innovative therapies with maximum benefits for patients.

### **About CEREP**

CEREP was founded by Thierry Jean in 1989. The company aims to optimise the drug discovery process.

CEREP offers the pharmaceuticals industry a range of solutions for cutting product development timescales and costs by identifying the molecules that are unsuitable for drug development at a very early stage. To achieve this, CEREP relies on a comprehensive technological platform focused on chemistry, biology and IT.

The company was listed on the *Nouveau Marché* of the Paris Stock Exchange on 18 February 1998 (SICOVAM code: 6179).

Contact:

Press Office  
CICOMMUNICATION  
Catherine Isnard  
Tel: +33-1 47 23 90 48  
E-mail: [cicom@cicommunication.fr](mailto:cicom@cicommunication.fr)