

**THE DEPARTMENT OF HEALTH OF THE GOVERNMENT  
OF THE UNITED KINGDOM**

**GOOD LABORATORY PRACTICE**

**STATEMENT OF COMPLIANCE  
IN ACCORDANCE WITH DIRECTIVE 2004/9/EC**

**TEST FACILITY**

CONCEPT LIFE SCIENCES INTERGRATED DISCOVERY  
AND DEVELOPMENT SERVICES LIMITED  
DISCOVERY PARK HOUSE  
RAMSGATE ROAD  
SANDWICH  
CT13 9ND  
UNITED KINGDOM

**TEST TYPE(S)**

Analytical/Clinical Chemistry

**DATE OF INSPECTION:** 06/10/2020

**DATE OF ISSUE:** 24/12/2020

An Inspection for compliance with the Principles of Good Laboratory Practice was carried out at the above named test facility as part of the UK Good Laboratory Practice Compliance Monitoring Programme.

This statement confirms that, on the date of issue, the UK Good Laboratory Practice Monitoring Authority were satisfied that the above named test facility was operating in compliance with the OECD Principles of Good Laboratory Practice.

This statement constitutes a Good Laboratory Practice Instrument (as defined in the UK Good Laboratory Practice Regulations 1999).

Issued by  
Mr Stephen Vinter  
Head, UK GLP Monitoring Authority

