



Department
of Health &
Social Care

**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 - DUNDEE
2 JAMES LINDSAY PLACE
DUNDEE TECHNOPOLE
DUNDEE
DD1 5JJ
UNITED KINGDOM

TEST TYPE(S)

Analytical Chemistry

DATE OF INSPECTION: 06/07/2020 - 21/07/2020

DATE OF ISSUE: 22/12/2020

Due to COVID-19 travel restrictions, a remote 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



Medicines & Healthcare products
Regulatory Agency

