



Department
of Health &
Social Care

THE DEPARTMENT OF HEALTH OF THE GOVERNMENT OF THE UNITED KINGDOM

GOOD LABORATORY PRACTICE

STATEMENT OF COMPLIANCE

TEST FACILITY

CONCEPT LIFE SCIENCES INTEGRATED DISCOVERY &
DEVELOPMENT SERVICES LIMITED
UNIT 69
LISTERHILLS SCIENCE PARK
CAMPUS ROAD
BRADFORD
BD7 1HR
UNITED KINGDOM

TEST TYPE(S)

Analytical/Clinical Chemistry
Physical/Chemical Testing

DATE OF INSPECTION: 10/02/2021

DATE OF ISSUE: 15/04/2021

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 UK Good Laboratory Practice Regulations, Statutory Instrument 1999 No. 3106 (as amended) which incorporate 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

