



Department
of Health

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

ASTERAND UK ACQUISITION LIMITED
2 ORCHARD ROAD
ROYSTON
SG8 5HD
UNITED KINGDOM

TEST TYPE(S)

Environmental Fate
Ecosystems
Other
In Vitro Safety Testing

DATE OF INSPECTION: 16/05/2017

DATE OF ISSUE: 30/10/2017

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
Dr Andrew J Gray
Head, UK GLP Monitoring Authority



Medicines & Healthcare products
Regulatory Agency



MHRA
Regulating Medicines and Medical Devices



Department
of Health

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

ASTERAND UK ACQUISITION LIMITED
2 ORCHARD ROAD
ROYSTON
SG8 5HD
UNITED KINGDOM

TEST TYPE(S)

Environmental Fate
Ecosystems
Other
In Vitro Safety Testing

DATE OF INSPECTION: 16/05/2017

DATE OF ISSUE: 30/10/2017

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
Dr Andrew J Gray
Head, UK GLP Monitoring Authority



Medicines & Healthcare products
Regulatory Agency



MHRA
Regulating Medicines and Medical Devices