



NSAI

Certificate of Registration of Quality Management System to ISO 9001:2015

The National Standards Authority of Ireland Inc. certifies that:

Quanterix
900 Middlesex Turnpike
Billerica, MA 01821
USA

has been assessed and deemed to comply with the requirements of the above standard in respect of the scope of operations given below:

Development, Manufacturing, Instrument and software Engineering and Engineering of assay instruments

Geraldine Larkin
Chief Executive Officer

Lisa Greenleaf
Technical Operations Officer

Registration Number: 19.8225
Certification Granted: May 08, 2019
Effective Date: November 07, 2022
Expiry Date: May 07, 2025



National Standards Authority of Ireland Inc., 20 Trafalgar Square, Nashua, New Hampshire, NH 03063, USA T +1 603 882 4412