



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

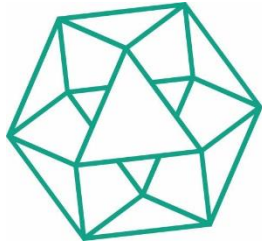
Geraldine Larkin
Chief Executive Officer

Lisa Greenleaf
Technical Operations Officer

Registration Number: 19.8225
Certification Granted: November 07, 2022
Effective Date: May 08, 2025
Expiry Date: May 07, 2028



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



NSAI

Certificate of Registration of Quality Management System to ISO 13485:2016

The National Standards Authority of Ireland certifies that:

Quanterix Corp.
900 Middlesex Turnpike,
Bldg.#1
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:

The design and development, manufacture, installation and servicing of reagents, test kits, instruments and software for neurological immunoassays.

Approved by:
Kevin Mullaney
Director of Certification

Registration Number: CM19.8685
Certification Granted: January 31, 2025
Effective Date: January 31, 2025
Expiry Date: January 30, 2028



National Standards Authority of Ireland, 1 Swift Square, Northwood, Santry, Dublin 9, Ireland T +353 1 807 3800
National Standards Authority of Ireland, 20 Trafalgar Square, Nashua, New Hampshire, NH 03063, USA T +1 603 882 4412