



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