



CERTIFICATE OF ASSURANCE

Revvity (UK) Ltd.

Llantrisant Business Park Llantrisant CF72 8YW

COMPLIES WITH THE REQUIREMENTS OF THE CYBER ESSENTIALS PLUS SCHEME

NAME OF ASSESSOR : Darren O'Regan

CERTIFICATE NUMBER : 502f22f2-d18d-499d-9f78-31debee756ad

DATE OF CERTIFICATION : 2024-05-01

PROFILE VERSION : 3.1 (Montpellier)

RECERTIFICATION DUE : 2025-05-01

SCOPE : Revvity (UK) Ltd. Clinical Laboratory Network excluding all other networks



SCAN QR CODE TO VERIFY THE AUTHENTICITY OF THIS CERTIFICATE

CERTIFICATION MARK



CERTIFICATION BODY



CYBER ESSENTIALS PARTNER



The Certificate certifies that the organisation was assessed as meeting the Cyber Essentials Plus implementation profile and thus that, at the time of testing, the organisations ICT defences were assessed as satisfactory against commodity based cyber attack. However, this Certificate does not in any way guarantee that the organisations defences will remain satisfactory against a cyber attack.