



COLLEGE of AMERICAN  
PATHOLOGISTS

## CERTIFICATE OF ACCREDITATION

**Revvity (UK) Ltd**  
**Revvity Omics**  
**Macclesfield, United Kingdom**  
**Jorune Balciuniene, PhD,ABMGG**

CAP#: 9602320

The organization named above meets all applicable standards for accreditation and is hereby accredited by the College of American Pathologists' Laboratory Accreditation Program. Reinspection should occur prior to **March 31, 2027** to maintain accreditation.

Accreditation does not automatically survive a change in director, ownership, or location and assumes that all interim requirements are met.

Kathleen G. Beavis, MD  
Chair, Accreditation Committee

Donald S. Karcher, MD, FCAP  
President, College of American Pathologists

