



COLLEGE of AMERICAN
PATHOLOGISTS

CERTIFICATE OF ISO 15189 ACCREDITATION

**Revvity (UK) Ltd. – Revvity Omics
Macclesfield, United Kingdom**

CAP # 9602320

The organization named above is granted this accreditation in accordance with the recognized International Standard ISO 15189:2012, Medical Laboratories – Requirements for quality and competence. This accreditation demonstrates competence for a defined scope and the operation of a laboratory quality management system.

Effective From August 4, 2023 through August 4, 2026

Vipul Trivedi, MD, FCAP
Chair, CAP 15189 Committee

Emily E. Volk, MD, FCAP
President, College of American Pathologists

The scope of this accreditation includes the Quality Management System and the discipline of Molecular Pathology.

