



COLLEGE of AMERICAN  
PATHOLOGISTS

# CERTIFICATE OF ACCREDITATION

**Revvity Omics  
Laboratory  
Pittsburgh, Pennsylvania  
Taraka R. Donti, PhD, MSc, BS, ABMGG**

CAP Number: 1286403  
AU-ID: 1178174  
CLIA Number: 39D0673919

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 **June 15, 2024** 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, Accreditation Committee Chair

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

