



COLLEGE of AMERICAN
PATHOLOGISTS

CERTIFICATE OF ACCREDITATION

Revvity Omics Inc
Laboratory
Pittsburgh, Pennsylvania
Taraka R. Donti, PhD,ABMGG

CAP#: 1286403
CLIA#: 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, 2028** to maintain accreditation.

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

Richard C. Friedberg, MD, PhD
Chair, Accreditation Committee

Qihui "Jim" Zhai, MD, FCAP
President, College of American Pathologists

