

**CENTERS FOR MEDICARE & MEDICAID SERVICES
CLINICAL LABORATORY IMPROVEMENT AMENDMENTS**

CERTIFICATE OF ACCREDITATION

LABORATORY NAME AND ADDRESS

REVVITY OMICS, INC
250 INDUSTRY DRIVE, SUITE 400
PITTSBURGH, PA 15275-1017

CLIA ID NUMBER

39D0673919

EFFECTIVE DATE

02/02/2026

LABORATORY DIRECTOR

DR. TARAKA R. DONTI

EXPIRATION DATE

02/01/2028

Pursuant to Section 353 of the Public Health Services Act (42 U.S.C. 263a) as revised by the Clinical Laboratory Improvement Amendments (CLIA), the above named laboratory located at the address shown hereon (and other approved locations) may accept human specimens for the purposes of performing laboratory examinations or procedures.

This certificate shall be valid until the expiration date above, but is subject to revocation, suspension, limitation, or other sanctions for violation of the Act or the regulations promulgated thereunder.



Gregg Brandush, Director
Division of Clinical Laboratory Improvement & Quality
Quality & Safety Oversight Group
Center for Clinical Standards and Quality

If you currently hold a Certificate of Compliance or Certificate of Accreditation, below is a list of the laboratory specialties/subspecialties you are certified to perform and their effective date:

LAB CERTIFICATION (CODE)	EFFECTIVE DATE	LAB CERTIFICATION (CODE)	EFFECTIVE DATE
MICROBIOLOGY - BACTERIOLOGY (110)	02/18/2014		
MICROBIOLOGY - VIROLOGY (140)	06/18/2010		
DIAGNOSTIC IMMUNOLOGY - GENERAL IMMUNOLOGY (220)	12/07/2025		
CHEMISTRY - ROUTINE CHEMISTRY (310)	03/29/2003		
CHEMISTRY - ENDOCRINOLOGY (330)	04/16/2008		
IMMUNOHEMATOLOGY - ANTIBODY DETECTION (NON-TRANSFUSION) (530)	12/12/2025		

**PLEASE CONTACT YOUR STATE AGENCY FOR ANY CHANGES TO YOUR CURRENT CERTIFICATE.
FOR MORE INFORMATION ABOUT CLIA, VISIT OUR WEBSITE AT WWW.CMS.GOV/CLIA.**