

CLINICAL LABORATORY PERMIT



pennsylvania
DEPARTMENT OF HEALTH

Pursuant to the act of September 26, 1951, P.L. 1539 as amended, a Permit to operate a Clinical Laboratory is hereby granted to:

Laboratory Identification Number: 22936A

AUTHORIZED CATEGORIES/TESTS:

Name and Director of Laboratory:

**CLINICAL CHEMISTRY
NON-SYPHILIS SEROLOGY
VIROLOGY**

**REVVITY OMICS, INC.
TARAKA R DONTI, PH.D.
250 INDUSTRY DRIVE, SUITE 400
PITTSBURGH, PA 15275-1017**

Owner:

REVVITY, INC

ISSUE DATE: August 15, 2024

DATE EXPIRES: August 15, 2025

Debra L. Bogen MD

**Debra L. Bogen, MD, FAAP
Acting Secretary of Health**

DISPLAY THIS CERTIFICATE PROMINENTLY

This permit is subject to revocation, suspension, or limitation for violation of the Act or the Regulations promulgated thereunder.

**REVVITY OMICS, INC.
TARAKA R DONTI, PH.D.
PO BOX 219
BRIDGEVILLE, PA 15017**