



MARYLAND DEPARTMENT OF HEALTH  
**Office of Health Care Quality**

*7120 Samuel Morse Drive, Second Floor, Columbia, MD 21046*

**NON-EXPIRING CLINICAL LABORATORY PERMIT**

**Type of Provider: Clinical Laboratory**

**License Number: 1072**

**Effective Date: 03/05/2024**

**Issued to: REVVITY OMICS INC  
250 INDUSTRY DRIVE, SUITE 400  
PITTSBURGH, PA 15275**

**Laboratory Director: Dr TARAKA DONTI**

**Owner of Laboratory: REVVITY INC.**

**Medical Laboratory Tests in the following disciplines may be performed:**

**Immunology:**

General Immunology

**Chemistry:**

Endocrinology, Routine

**Molecular Biology:**

Nucleic Acid Probes

**Genetics:**

Molecular

This license is issued by the Maryland Department of Health in accordance with the applicable provisions of Maryland Code Ann., Health General Article and its rules and regulations. This license is not transferable.

*Patricia Tomsko Nay, MD*

Patricia Tomsko Nay, MD

Executive Director

Office of Health Care Quality

Falsification of a license shall subject the perpetrator to criminal prosecution and the imposition of civil fines.