

# New York State Department of Health

PFI: 4733

## Clinical Laboratory Permit

CLIA: 39D0673919

Revvity Omics Inc

250 Industry Drive Suite 400

Pittsburgh PA 15275-1017

Director:  
Taraka R Donti, Ph.D.

Owner:  
Revvity Omics, Inc

is hereby authorized to perform laboratory procedures at the above location in the following categories in accordance with Article 5, Title V, Section 575 of the Public Health Law. This permit shall become void upon a change in the director, owner or location of the laboratory, and an application for a new permit shall be made to the Department.

*Diagnostic Immunology*  
*Diagnostic Services Serology*  
*Endocrinology*

*Genetic Testing*  
*Biochemistry*  
*(limited to screening and monitoring tests only)*  
*Molecular*

*Virology*  
*(limited to molecular methods)*

Amended

Effective Date: August 1, 2023

Expiration Date: June 30, 2024

Subject to Revocation

Permit Not Transferable

POST CONSPICUOUSLY

Serial: LAP 183140