

New York State Department of Health

PFL: 9325

Clinical Laboratory Permit

CLIA: 22DD2093470

Exosome Diagnostics, Inc.

266 Second Ave, Suite 200

Waltham MA 02451

Director:
Jordan S. Laser, M.D.

Owner:
Exosome Diagnostics, 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.

*Oncology
Molecular and Cellular Tumor Markers*

Renewal

Effective Date: July 1, 2024

Expiration Date: June 30, 2025

Subject to Revocation
Permit Not Transferable

POST CONSPICUOUSLY

Serial: LAP 189419