



State of Rhode Island
DEPARTMENT OF HEALTH
Center for Health Facilities Regulation

This is to certify that REVVITY OMICS, INC
250 INDUSTRY DRIVE SUITE 400 PITTSBURGH PA 15275
License Number: LCO01175

is hereby authorized to conduct and maintain an Out of State Clinical Laboratory in conformity with RIGL C23-16.2 and the standards, rules and regulations prescribed thereunder. This license is subject to biennial renewal unless sooner suspended or revoked for cause. The name on this license is the common name under which the licensee does business and may not reflect the legal license holder. Please call (401) 222-2566 for more information.

APPROVED SPECIALTY (IES)

MICROBIOLOGY, Virology, DIAGNOSTIC IMMUNOLOGY, General Immunology, CHEMISTRY, Routine Chemistry, Endocrinology, CLINICAL GENETICS, Molecular Genetics, Biochemical Genetics,

Expires: 12/30/2027

Issued: 01/19/2018

License Owner: REVVITY, INC- Corporation