

Revvity Omics, Inc.

Pittsburgh, PA

has been Accredited by




The Joint Commission


Which has surveyed this organization and found it to meet the requirements for the
Laboratory Accreditation Program

May 5, 2022

Accreditation is customarily valid for up to 24 months.


Jane Englebright, PhD, RN, CENP, FAAN
Chair, Board of Commissioners

ID #372985
Print/Reprint Date: 09/22/2023


Jonathan B. Perlin, MD, PhD, MSHA, MACP, FACMI
President and Chief Executive Officer

The Joint Commission is an independent, not-for-profit national body that oversees the safety and quality of health care and other services provided in accredited organizations. Information about accredited organizations may be provided directly to The Joint Commission at 1-800-994-6610. Information regarding accreditation and the accreditation performance of individual organizations can be obtained through The Joint Commission's web site at www.jointcommission.org.

