



Has Certified That  
Revvity Omics, Inc.

COLA ID #: 15860  
CLIA ID #: 39D0673919

Has met all criteria for  
**Laboratory Accreditation**

12/13/2023

This certificate expires two years from this date

A handwritten signature in black ink that reads 'Keith E. Davis MD'.

Chair of the Board

REV 0923

A handwritten signature in black ink that reads 'Nancy Stratton'.

Chief Executive Officer