

Product Number:	103360
Lot Number:	508412
Expiration:	25-Mar-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.96	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

Name

Title

Signature/Date

  27 MAR 2026

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

techsupport@quanterix.com

©Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures