

Product Number:	103360
Lot Number:	525904
Expiration:	16-Sep-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Kathleen Cornell

Quality Control
Supervisor

[Signature]
19 Sep 2025

Name

Title

Signature/Date

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