

Product Number:	102914
Lot Number:	523402
Expiration:	22-Aug-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.01	7.95 - 8.05	PASS

Review/Approval

Heidi Caswell
Name

Director, QC
Title

Heidi Caswell 08 Sep 2025
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