

Product Number:	101954
Lot Number:	523809
Expiration:	26-Feb-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.81	7.75 - 7.85	PASS

Review/Approval

Heidi Caswell

Director, QC

Heidi Caswell 08 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