

Product Number:	101357
Lot Number:	509413
Expiration:	4-Oct-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.18	6.18 - 6.22	PASS

Review/Approval

<i>Heidi Caswell</i>	<i>Director, QC</i>	<i>Heidi Caswell</i> 07Jul2025
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