

Product Number:	103158
Lot Number:	506220
Control Lot Number:	506220A
Expiration:	3-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.38	7.35 - 7.45	PASS

Review/Approval

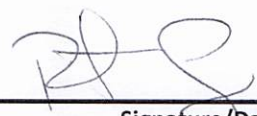
Robert Somogyi

QC Manager

Name

Title

Signature/Date

 18 MAR 2025

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