

Product Number:	101954
Lot Number:	414208
Expiration:	21-Nov-2025
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager

 28 MAY 2024

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