

Product Number:	101476
Lot Number:	508404
Expiration:	25-Sep-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.96	8.95 - 9.05	PASS

Review/Approval

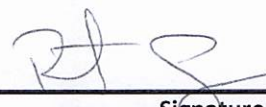
Robert Somogyi

QC Manager

Name

Title

Signature/Date

 27 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