

Product Number:	102914
Lot Number:	434018
Expiration:	5-Dec-2025
Storage:	2-8°C

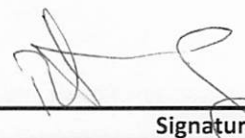
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.96	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title


Signature/Date

14 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