

Product Number:	102002
Lot Number:	506407
Expiration:	5-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

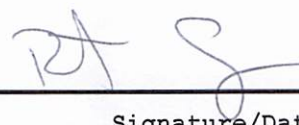
Robert Somogyi

QC Manager

Name

Title

Signature/Date

 26 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