

Product Number:	100458
Lot Number:	526507
Expiration:	22-Mar-2027
Storage:	2-8 °C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature] 26 Sep 2025

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