

Product Number:	100457
Lot Number:	521011
Expiration:	29-Jan-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

[Signature]
18 Aug 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