

Product Number:	101476
Lot Number:	431106
Expiration:	6-Nov-2025
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Heidi Caswell

Director, QC

H. Caswell 26 Nov 2024

Name

Title

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