

Product Number:	101353
Lot Number:	404108
Expiration:	10-Aug-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.41	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 05 MAR 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