

Product Number:	103361
Lot Number:	327910
Expiration:	6-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.02	7.95 - 8.05	PASS

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 11 OCT 2023
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