

Product Number:	103361
Lot Number:	403711
Expiration:	6-Feb-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.04	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 14 FEB 2024

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