

Product Number:	102907
Lot Number:	503942
Expiration:	19-Jul-2024
Storage:	2 - 8°C

Component	Part Number	Lot Number	Characteristic	Pass / Fail
Lysate Diluent A Reagent	102908	320006	pH meter	Pass
Lysate Diluent B Reagent	102910	320008		Pass
Lysate Diluent C Reagent	102912	320102		Pass
Lysate Diluent D Reagent	102914	322311		Pass

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan 29 AUG 2023

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