

Product Number:	101477
Lot Number:	503961
Expiration:	12-May-2024
Storage:	2-8°C

Component	Part Number	Lot Number	Control Number
Sample Diluent A	101472	315310	315310A
Sample Diluent B	101473	319914	319914A
Sample Diluent C	101474	317204	317204A
Sample Diluent D	101475	313205	313205A
Sample Diluent E ¹	101476	322906	n/a

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45
Characteristic	Result	pH Target
pH	Pass	8.95 - 9.05 ¹

¹ Sample Diluent E (101476) pH target differs from the all other components
Note: Result corresponds to each component

Review/Approval

Sonia Khillar
Name

QC Manager
Title

Sonia Khillar 06SEP2023
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