

Product Number:	101477
Lot Number:	503687
Expiration:	23-Apr-2023
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	212415
Sample Diluent B	101473	211303
Sample Diluent C	101474	214901
Sample Diluent D	101475	215507
Sample Diluent E ¹	101476	214903

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 Khillan
Name

QC Manager
Title

Sonia Khillan / 30 AUG 2022
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