

Product Number:	101477
Lot Number:	504020
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	328202	N/A
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 Khillan

Senior QC Analyst

Sonia Khillan 19 OCT 2023

Name

Title

Signature/Date