

Product Number:	101477
Lot Number:	503886
Expiration:	23-Feb-2024
Storage:	2-8°C

Component	Part Number	Lot Number	Control Number
Sample Diluent A	101472	306618	n/a
Sample Diluent B	101473	308218	n/a
Sample Diluent C	101474	312906	312906A
Sample Diluent D	101475	313204	313204A
Sample Diluent E ¹	101476	305404	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
Name

QC Manager
Title

Sonia 12/11- 30 MAY 2023
Signature/Date