

Product Number:	101477
Lot Number:	503776
Expiration:	29-May-2023
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	230509
Sample Diluent B	101473	230115
Sample Diluent C	101474	234812
Sample Diluent D	101475	233302
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 12/12/2022
Signature/Date