

Product Number:	101477
Lot Number:	503801
Expiration:	28-Oct-2023
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	230509
Sample Diluent B	101473	230115
Sample Diluent C	101474	301712
Sample Diluent D	101475	233302
Sample Diluent E ¹	101476	234905

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

Matt Koulouris
Name

QC Supervisor
Title

[Signature] 07Feb2023
Signature/Date