

Product Number:	101477
Lot Number:	503528
Expiration:	4-May-2023
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	212415
Sample Diluent B	101473	215103
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

Matt Smerit
Name

QC
Title

Mda 27 Jun 2022
Signature/Date