

Product Number:	101477
Lot Number:	503397
Expiration:	30-Jul-2022
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	122011
Sample Diluent B	101473	121118
Sample Diluent C	101474	204908
Sample Diluent D	101475	123210
Sample Diluent E ¹	101476	134401

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

ROBERT SOMOSYI AS
16 MAR 2022

Shivani Goel

QC Manager

Name

Title

Signature/Date
16 MAR 2022