

Product Number:	101477
Lot Number:	503426
Expiration:	8-Aug-2022
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	122011
Sample Diluent B	101473	133937
Sample Diluent C	101474	204908
Sample Diluent D	101475	135005
Sample Diluent E ¹	101476	201811

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

Shivani Goel

QC Manager

Name

Title

Shivani Goel 04Apr2022

Signature/Date