

Product Number:	101477
Lot Number:	503221
Expiration:	31-Mar-2022
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	109023
Sample Diluent B	101473	115231
Sample Diluent C	101474	124003
Sample Diluent D	101475	123210
Sample Diluent E	101476	123306

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Note: Result corresponds to each component

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 17 Nov 2021
Signature/Date