

Product Number:	101477
Lot Number:	503730
Expiration:	4-Jun-2023
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	230509
Sample Diluent B	101473	227215
Sample Diluent C	101474	227106
Sample Diluent D	101475	215507
Sample Diluent E ¹	101476	224211

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

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 08NOV2022
Signature/Date