

Product Number:	101477
Lot Number:	503991
Expiration:	18-Jul-2024
Storage:	2-8°C

Component	Part Number	Lot Number	Control Number
Sample Diluent A	101472	324510	N/A
Sample Diluent B	101473	319918	319918A
Sample Diluent C	101474	326408	N/A
Sample Diluent D	101475	326106	N/A
Sample Diluent E ¹	101476	326309	N/A

Laboratory Analysis

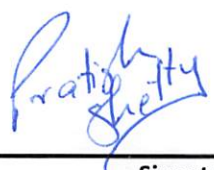
Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45
Characteristic	Result	pH Target
pH	Pass	8.95 - 9.05 ¹

¹ Sample Diluent E (101476) pH target differs from the all other components

Note: Result corresponds to each component

Review/Approval





Name Title Signature/Date

Pratishk Shetty SUPERVISOR 27 SEP 2023