

Product Number:	101477
Lot Number:	504175
Expiration:	14-Feb-2025
Storage:	2-8°C


Component	Part Number	Lot Number
Sample Diluent A	101472	402615
Sample Diluent B	101473	401614
Sample Diluent C	101474	404514
Sample Diluent D	101475	404416
Sample Diluent E ¹	101476	404609

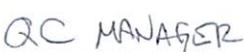
Laboratory Analysis


Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45
Characteristic	Result	pH Target
pH Measurement	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