

Product Number:	101477
Lot Number:	504452
Expiration:	20-May-2025
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	407815
Sample Diluent B	101473	401622
Sample Diluent C	101474	429508
Sample Diluent D	101475	425610
Sample Diluent E ¹	101476	414115

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

TAM LEE DOC CONTROL II, QA *Tam Lee* 15 NOV 2024
 Name Title Signature/Date