

Product Number:	101477
Lot Number:	504716
Expiration:	16-Jun-2026
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	514708
Sample Diluent B	101473	503107
Sample Diluent C	101474	519805
Sample Diluent D	101475	521907
Sample Diluent E ¹	101476	516704

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

Dhwani Patel

QA Specialist II, QA

 03 Sep 2025

NameTitleSignature/Date