

Product Number:	101477
Lot Number:	504737
Expiration:	5-Aug-2026
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	514710
Sample Diluent B	101473	521616
Sample Diluent C	101474	521712
Sample Diluent D	101475	521907
Sample Diluent E ¹	101476	526506

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

SDR / 26 Sep 2025

Name

Title

Signature/Date