

Product Number:	101477
Lot Number:	504694
Expiration:	22-Apr-2026
Storage:	2-8°C

Component	Part Number	Lot Number	Control Lot Number
Sample Diluent A	101472	501527	501527A
Sample Diluent B	101473	503107	N/A
Sample Diluent C	101474	511201	N/A
Sample Diluent D	101475	521907	N/A
Sample Diluent E ¹	101476	516704	N/A

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

Dilish Patel
Name

Sr. QA Specialist
Title

21 *19 AUG 2025*
Signature/Date