

Product Number:	101477
Lot Number:	504605
Expiration:	22-Apr-2026
Storage:	2-8°C

Component	Part Number	Lot Number	Control Lot Number
Sample Diluent A	101472	501524	501524A
Sample Diluent B	101473	506221	N/A
Sample Diluent C	101474	511201	N/A
Sample Diluent D	101475	510607	N/A
Sample Diluent E ¹	101476	508404	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

Dipesh Patel

Sec. Doc. Control

29 APR 2025

Name

Title

Signature/Date