

Product Number:	101477
Lot Number:	504572
Expiration:	6-Nov-2025
Storage:	2-8°C

Component	Part Number	Lot Number	Control Lot Number
Sample Diluent A	101472	501524	501524A
Sample Diluent B	101473	426818	N/A
Sample Diluent C	101474	505208	N/A
Sample Diluent D	101475	425610	N/A
Sample Diluent E ¹	101476	431106	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

Name

Sen. Doc control - QA

Title

NT 102 APR 2025

Signature/Date