

Product Number:	101477
Lot Number:	504160
Expiration:	13-Oct-2024
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	402615
Sample Diluent B	101473	401614
Sample Diluent C	101474	404514
Sample Diluent D	101475	404416
Sample Diluent E ¹	101476	328612

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

ROBERT SOMOGYI

QC MANAGER

[Signature] 29 FEB 2024

Name

Title

Signature/Date