

Product Number:	101477
Lot Number:	504087
Expiration:	23-Mar-2024
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	329410
Sample Diluent B	101473	308223
Sample Diluent C	101474	334013
Sample Diluent D	101475	326107
Sample Diluent E ¹	101476	326310

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

Sonia Khillan

Name

Senior QC Analyst

Title

Sonia 12.11.23 13 DEC 2023

Signature/Date