

Product Number:	101477
Lot Number:	504488
Expiration:	16-Jul-2025
Storage:	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	415509
Sample Diluent B	101473	401614
Sample Diluent C	101474	434504
Sample Diluent D	101475	425610
Sample Diluent E <sup>1</sup>	101476	431106

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45
Characteristic	Result	pH Target
pH Measurement	Pass	8.95 - 9.05 <sup>1</sup>

<sup>1</sup> Sample Diluent E (101476) pH target differs from the all other components

Note: Result corresponds to each component

Review/Approval

Dipesh Patel  
Name

Doc control TE QA  
Title

CT 27 DEC 2024  
Signature/Date