

<b>Product Number:</b>	101477
<b>Lot Number:</b>	503757
<b>Expiration:</b>	29-Aug-2023
<b>Storage:</b>	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	230509
Sample Diluent B	101473	227215
Sample Diluent C	101474	227106
Sample Diluent D	101475	233302
Sample Diluent E <sup>1</sup>	101476	224211

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45
Characteristic	Result	pH Target
pH	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

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan 08DEC2022*  
Signature/Date