

<b>Product Number:</b>	101477
<b>Lot Number:</b>	502952
<b>Expiration:</b>	15-Feb-2022
<b>Storage:</b>	2-8°C

Component	Part Number	Lot Number
Sample Diluent A	101472	109023
Sample Diluent B	101473	104617
Sample Diluent C	101474	109211
Sample Diluent D	101475	107808
Sample Diluent E	101476	109213

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Note: Result corresponds to each component

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 01Jun2021  
Signature/Date