

<b>Product Number:</b>	103847
<b>Lot Number:</b>	104917
<b>Expiration:</b>	18-Feb-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.35-7.45

Review/Approval

*Shivani Goel*  
Name

*QC Manager*  
Title

*Shivani Goel 12-Mar-2021*  
Signature/Date