

<b>Product Number:</b>	103847
<b>Lot Number:</b>	206810
<b>Expiration:</b>	9-Mar-2023
<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

QC Manager

*Shivani Goel 04Apr 2022*

Name

Title

Signature/Date