

<b>Product Number:</b>	101631
<b>Lot Number:</b>	116810
<b>Expiration:</b>	17-Jun-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.30-7.50

Review/Approval

*Shivani Goel*

Name

QC

Title

*Shivani Goel 07/01/2021*

Signature/Date