

<b>Product Number:</b>	101631
<b>Lot Number:</b>	034205
<b>Expiration:</b>	7-Dec-2021
<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

Junpil Park

Supervisor, QC

 15 Feb 2021

Name

Title

Signature/Date