

<b>Product Number:</b>	103847
<b>Lot Number:</b>	029315
<b>Expiration:</b>	19-Oct-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  
Name

QC Supervisor  
Title

 14Dec2020  
Signature/Date