

<b>Product Number:</b>	103847
<b>Lot Number:</b>	117119
<b>Expiration:</b>	20-Jun-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  
Title

Shivani Goel 29 Jul 2021  
Signature/Date