

<b>Product Number:</b>	102252
<b>Lot Number:</b>	123909
<b>Expiration:</b>	27-Aug-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 24Feb2022  
Signature/Date