

<b>Product Number:</b>	102910
<b>Lot Number:</b>	127408
<b>Expiration:</b>	1-Oct-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.90-8.10

Review/Approval

Shivam Goel  
Name

QC  
Title

Shivam Goel 12Oct2021  
Signature/Date