

<b>Product Number:</b>	102104
<b>Lot Number:</b>	110015
<b>Expiration:</b>	10-Apr-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 21Jan2022*

Signature/Date