

<b>Product Number:</b>	102912
<b>Lot Number:</b>	235008
<b>Expiration:</b>	16-Dec-2023
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.95 - 8.05

Review/Approval

*Sonia Khillan*

Name

*QC Manager*

Title

*Sonia Khillan 103 JAN 2023*

Signature/Date