

<b>Product Number:</b>	101362
<b>Lot Number:</b>	300612
<b>Expiration:</b>	6-Jan-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	7.75 - 7.85

Review/Approval

*Sonia Khillan*

*QC Manager*

*Sonia Khillan 01MAY2023*

Name

Title

Signature/Date