

<b>Product Number:</b>	100458
<b>Lot Number:</b>	301313
<b>Expiration:</b>	13-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*

Name

*QC manager*

Title

*Sonia Khillan / 31 JAN 2023*

Signature/Date