

<b>Product Number:</b>	100458
<b>Lot Number:</b>	232109
<b>Expiration:</b>	17-Nov-2023
<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 Khilkin*

Name

*QC Manager*

Title

*Sonia Khilkin / 22 Nov 2022*

Signature/Date