

<b>Product Number:</b>	100458
<b>Lot Number:</b>	220106
<b>Expiration:</b>	20-Jul-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 Khillan*

Name

QC

Title

*Sonia Khillan / 27 Jul 2022*

Signature/Date