

<b>Product Number:</b>	103359
<b>Lot Number:</b>	223613
<b>Expiration:</b>	24-Aug-2023
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

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

Review/Approval

*Sonia Khillan*

Name

QC

Title

*Sonia Khillan / 29 AUG 2022*

Signature/Date