

<b>Product Number:</b>	101357
<b>Lot Number:</b>	223710
<b>Expiration:</b>	25-Aug-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	6.18 - 6.22

Review/Approval

*Sonia Khillan*

Name

QC

Title

*Sonia Khillan 129 AUG 2022*

Signature/Date