

<b>Product Number:</b>	103847
<b>Lot Number:</b>	322711
<b>Expiration:</b>	15-Aug-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH	7.36	7.35 - 7.45	PASS

Review/Approval

*Sonia Killian*

Name

*QC Manager*

Title

*Sonia Killian 26SEP2023*

Signature/Date