

<b>Product Number:</b>	103847
<b>Lot Number:</b>	226216
<b>Expiration:</b>	19-Sep-2023
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan QC Manager Sonia 12/16/25 OCT 2022

Name

Title

Signature/Date