

<b>Product Number:</b>	101473
<b>Lot Number:</b>	227215
<b>Expiration:</b>	29-Aug-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

<u>Sonia Khillan</u>	<u>QC Manager</u>	<u></u>	<u>06SEP2022</u>
Name	Title	Signature/Date	