

<b>Product Number:</b>	103172
<b>Lot Number:</b>	231316
<b>Expiration:</b>	9-Nov-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

*Matt Koulouris*

Name

QC Supervisor

Title

*[Signature]* 15-Feb-2023

Signature/Date