

<b>Product Number:</b>	101472
<b>Lot Number:</b>	234802
<b>Expiration:</b>	14-Dec-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan / 24 JAN 2023

Signature/Date