

<b>Product Number:</b>	101355
<b>Lot Number:</b>	225111
<b>Expiration:</b>	8-Sep-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 12/16 / 13 SEP 2022  
Signature/Date