

<b>Product Number:</b>	101956
<b>Lot Number:</b>	301312
<b>Expiration:</b>	13-Jan-2024
<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

*[Signature] 12/16/31 JAN 2023*

Signature/Date