

<b>Product Number:</b>	101954
<b>Lot Number:</b>	327613
<b>Expiration:</b>	3-Oct-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH Measurement	7.75	7.75 - 7.85	PASS

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan 11 OCT 2023*  
Signature/Date