

<b>Product Number:</b>	101954
<b>Lot Number:</b>	318007
<b>Control Number:</b>	318007A
<b>Expiration:</b>	29-Jun-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval      Sonia Khillan      QC Manager      Sonia Khillan 12 JUL 2023

Name                                  Title                                  Signature/Date