

<b>Product Number:</b>	100458
<b>Lot Number:</b>	318009
<b>Control Number:</b>	318009A
<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 19 JUL 2023  
Name Title Signature/Date