

<b>Product Number:</b>	100458
<b>Lot Number:</b>	235305
<b>Expiration:</b>	19-Dec-2023
<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  
Name

QC Manager  
Title

Sonia Khillan / 12/23/2022  
Signature/Date