

<b>Product Number:</b>	103158
<b>Lot Number:</b>	301311
<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 Khillsu*  
Name

*QC Manager*  
Title

*Sonia R. / 13 JAN 2023*  
Signature/Date