

<b>Product Number:</b>	100453
<b>Lot Number:</b>	303210
<b>Expiration:</b>	1-Feb-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	6.18 - 6.22

Review/Approval

*Sonia Khillan*  
Name

QC

Title

*Sonia Khillan / 13 FEB 2023*  
Signature/Date