

<b>Product Number:</b>	101353
<b>Lot Number:</b>	217217
<b>Expiration:</b>	21-Jun-2023
<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 Khillan  
Name

QC Manager  
Title

Sonia 12/11/2022  
Signature/Date