

<b>Product Number:</b>	101954
<b>Lot Number:</b>	208805
<b>Expiration:</b>	29-Mar-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

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 01 Apr 2022  
Signature/Date