

<b>Product Number:</b>	101352
<b>Lot Number:</b>	111608
<b>Expiration:</b>	26-Apr-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
pH Measurement	PASS

Review/Approval

*Shivani Gael*  
Name

*QC Manager*  
Title

*Shivani Gael 03.May 2021*  
Signature/Date