

<b>Product Number:</b>	101352
<b>Lot Number:</b>	117225
<b>Expiration:</b>	21-Jun-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 04 Aug 2021  
Signature/Date