

<b>Product Number:</b>	101353
<b>Lot Number:</b>	131516
<b>Expiration:</b>	11-Nov-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

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

Review/Approval

Shivani Goel  
Name

Quality Control  
Title

Shivani Goel 08 Dec 2021  
Signature/Date