

<b>Product Number:</b>	101357
<b>Lot Number:</b>	134608
<b>Expiration:</b>	12-Dec-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 20Dec2021*  
Signature/Date