

<b>Product Number:</b>	101355
<b>Lot Number:</b>	122510
<b>Expiration:</b>	13-Aug-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval Shivani Goel QC Shivani Goel 23 Aug 2021  
Name Title Signature/Date