

<b>Product Number:</b>	101352
<b>Lot Number:</b>	031122
<b>Expiration:</b>	6-Nov-2021
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

**Review/Approval**

Shivani Goel

Manager, QC

*Shivani Goel 10Feb2021*

**Name**

**Title**

**Signature/Date**