

<b>Product Number:</b>	101362
<b>Lot Number:</b>	026617
<b>Expiration:</b>	22-Sep-2021
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

**Review/Approval**

Shivani Goel

QC Manager

**Name**

**Title**

*Shivani Goel (6 Oct 2020)*  
**Signature/Date**