

<b>Product Number:</b>	103158
<b>Lot Number:</b>	029013
<b>Expiration:</b>	16-Oct-2021
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Shivani Goel

Manager, QC

Name

Title

Shivani Goel 24 Nov 2020

Signature/Date