

<b>Product Number:</b>	100458
<b>Lot Number:</b>	110117
<b>Expiration:</b>	11-Apr-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

*Shivani Goel*

Name

*QC Manager*

Title

*Shivani Goel 03 May 2021*

Signature/Date