

<b>Product Number:</b>	101362
<b>Lot Number:</b>	110119
<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

QC Manager

Shivani Goel 24 May 2021

Name

Title

Signature/Date