

<b>Product Number:</b>	100453
<b>Lot Number:</b>	100706
<b>Expiration:</b>	7-Jan-2022
<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 10 Feb 2021*

Name

Title

Signature/Date