

<b>Product Number:</b>	100453
<b>Lot Number:</b>	127413
<b>Expiration:</b>	1-Oct-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

*Shivani Goel*  
Name

QC

Title

*Shivani Goel 21-Jan-2022*  
Signature/Date