

<b>Product Number:</b>	100376
<b>Lot Number:</b>	203113
<b>Expiration:</b>	31-Jan-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Shivani Goel  
Name

QC

Title

Shwami Goel 15Feb2022  
Signature/Date