

<b>Product Number:</b>	100376
<b>Lot Number:</b>	211404
<b>Expiration:</b>	24-Apr-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Richa Shrestha

QC Supervisor

Name

Title

 26Apr2022  
Signature/Date