

<b>Product Number:</b>	101362
<b>Lot Number:</b>	101605
<b>Expiration:</b>	16-Jan-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Junpil Park

Supervisor, QC



22 Feb 2021

Name

Title

Signature/Date