

<b>Product Number:</b>	101356
<b>Lot Number:</b>	024109
<b>Expiration:</b>	28-Aug-2021
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Junpil Park

Supervisor, QC

 30 Dec 2020

Name

Title

Signature/Date