

<b>Product Number:</b>	101955
<b>Lot Number:</b>	104721
<b>Expiration:</b>	16-Feb-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

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

Review/Approval

Junpil Park

Quality Control, Supervisor

 04Mar2021

Name

Title

Signature/Date