

<b>Product Number:</b>	101472
<b>Lot Number:</b>	117712
<b>Expiration:</b>	26-Jun-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.20-7.60

Review/Approval

*Junpil Park*

Name

*QC Supervisor*

Title

*[Signature]* *01 Sep 2021*

Signature/Date