

<b>Product Number:</b>	101472
<b>Lot Number:</b>	109023
<b>Expiration:</b>	31-Mar-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.35-7.45

Review/Approval

*Jungil Park*

Name

QC Supervisor

Title

*[Signature]* 22 Apr 2021

Signature/Date