

<b>Product Number:</b>	103359
<b>Lot Number:</b>	124001
<b>Expiration:</b>	28-Aug-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.90-8.10

Review/Approval

*Junpil Park*

Name

QC Supervisor

Title

*[Signature]* 01sep2021

Signature/Date