

<b>Product Number:</b>	101556
<b>Lot Number:</b>	031524
<b>Expiration:</b>	10-Nov-2021
<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

Junpil Park  
Name

QC Supervisor  
Title

 30 Dec 2020  
Signature/Date