

<b>Product Number:</b>	102912
<b>Lot Number:</b>	102509
<b>Expiration:</b>	25-Jan-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.95-8.05

Review/Approval

Junpil Park  
Name

Supervisor, Quality Control  
Title

 02 Feb 2021  
Signature/Date