

<b>Product Number:</b>	101356
<b>Lot Number:</b>	110521
<b>Expiration:</b>	15-Apr-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
pH Measurement	PASS

Review/Approval

*J. Park*  
Name

QC SUPERVISOR  
Title

*[Signature]* 19 Apr 2021  
Signature/Date