

<b>Product Number:</b>	101357
<b>Lot Number:</b>	216704
<b>Expiration:</b>	16-Jun-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	6.18 - 6.22

Review/Approval

*Matt Smart*

Name

QC

Title

*MAT 22 Jun 2022*

Signature/Date