

<b>Product Number:</b>	102002
<b>Lot Number:</b>	225606
<b>Expiration:</b>	13-Sep-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

*Matt Smart*

*QC*

*MSA 19 Sep 2022*

Name

Title

Signature/Date