

Product Number:	103847
Lot Number:	216002
Expiration:	9-Jun-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

*Matt Smart*

Name

QC

Title

*MSA* 13Jun2022

Signature/Date