

Product Number:	101475
Lot Number:	215507
Expiration:	4-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