

Product Number:	103359
Lot Number:	217311
Expiration:	22-Jun-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

*Matt Smart*  
Name

QC

Title

*[Signature]* 27 Jun 2022  
Signature/Date