

Product Number:	102908
Lot Number:	215504
Expiration:	4-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

MA 14 Jun 2022  
Signature/Date