

Product Number:	103360
Lot Number:	210920
Expiration:	19-Apr-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.90 - 8.10

Review/Approval

Matt Smart  
Name

QC  
Title

MJA 19 May 2022  
Signature/Date