

Product Number:	103358
Lot Number:	210905
Expiration:	19-Apr-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel 20Apr2022*  
Signature/Date