

Product Number:	102908
Lot Number:	210904
Expiration:	19-Apr-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 20 Apr

Name

Title

Signature/Date

2022