

Product Number:	103358
Lot Number:	205703
Expiration:	26-Feb-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 07Mar2022*

Name

Title

Signature/Date