

Product Number:	103361
Lot Number:	205712
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 07 Mar 2022*

Name

Title

Signature/Date