

Product Number:	102912
Lot Number:	205708
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