

Product Number:	103359
Lot Number:	210313
Expiration:	13-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 2022

Name

Title

Signature/Date