

Product Number:	101476
Lot Number:	210308
Expiration:	13-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 14Apr2022

Name

Title

Signature/Date