

Product Number:	101358
Lot Number:	209003
Expiration:	31-Mar-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 04Apr2022

Name

Title

Signature/Date