

Product Number:	103358
Lot Number:	213308
Expiration:	13-May-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 17 May 2022

Name

Title

Signature/Date