

Product Number:	103358
Lot Number:	109402
Expiration:	4-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel

Name

Quality Control

Title

Shivani Goel 14Apr2021

Signature/Date