

Product Number:	103359
Lot Number:	211308
Expiration:	23-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Richa Shrestha

Name

QC Supervisor

Title

[Signature] 26Apr2022

Signature/Date