

Product Number:	101473
Lot Number:	211303
Expiration:	23-Apr-2023
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Richa Shrestha

QC Supervisor

 27 Apr 2022

Name

Title

Signature/Date