

Product Number:	101358
Lot Number:	203014
Expiration:	30-Jan-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval


Richa Shrestha

QC Supervisor

Name

Title

Signature/Date

 21A382022