

Product Number:	103847
Lot Number:	110803
Expiration:	18-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park

Name

QC Supervisor

Title

[Signature] 22 Apr 2024

Signature/Date