

Product Number:	102620
Lot Number:	121310
Expiration:	1-Aug-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Jumpil Park

Name

QC Supervisor

Title

[Signature] 01sep2021

Signature/Date