

Product Number:	101475
Lot Number:	123210
Expiration:	20-Aug-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]

Signature/Date

01 Sep 2021