

Product Number:	101474
Lot Number:	109211
Expiration:	2-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Tunpil Park

Name

QC

Supervisor

Title

[Signature] 22 Apr 2021

Signature/Date