

Product Number:	103704
Lot Number:	101207
Expiration:	12-Jul-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park

Name

QC Supervisor

Title

 18 Feb 2021

Signature/Date