

Product Number:	101473
Lot Number:	104617
Expiration:	15-Feb-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



24 Feb 2021

Signature/Date