

Product Number:	101472
Lot Number:	033714
Expiration:	2-Dec-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park

QC Supervisor

 15 Feb 2021

Name

Title

Signature/Date