

Product Number:	100460
Lot Number:	218109
Expiration:	30-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

ROBERT SIMOIA

Name

QC

Title

 11/06/2022

Signature/Date