

Product Number:	100460
Lot Number:	215211
Expiration:	1-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan 08 JUL 2022

Signature/Date