

Product Number:	101475
Lot Number:	233302
Expiration:	29-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan / 07 DEC 2022

Name

Title

Signature/Date