

Product Number:	102908
Lot Number:	218909
Expiration:	8-Jul-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia Khillan / 25 JUL 2022
Signature/Date