

Product Number:	103358
Lot Number:	223409
Expiration:	22-Aug-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 / 24 AUG 2022
Signature/Date