

Product Number:	103361
Lot Number:	219212
Expiration:	11-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 / 13 JUL 2022
Signature/Date