

Product Number:	102286
Lot Number:	217603
Expiration:	25-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 / 25AUG2022
Signature/Date