

Product Number:	102914
Lot Number:	219211
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