

Product Number:	102914
Lot Number:	223615
Expiration:	24-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sania Khillan

Name

QC

Title

Sania Khillan / 29 AUG 2022

Signature/Date