

Product Number:	103847
Lot Number:	211315
Expiration:	23-Apr-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 / 03AVG2022

Signature/Date