

Product Number:	103847
Lot Number:	219601
Expiration:	15-Jul-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

QC

Sonia Khillan 12/21 JUL 2022

Name

Title

Signature/Date