

Product Number:	103847
Lot Number:	211813
Expiration:	28-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Matt Smart

QC

MMA 15 Aug 2022

Name

Title

Signature/Date