

Product Number:	103359
Lot Number:	214702
Expiration:	27-May-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Matt Smart

Name

QC

Title

M A 31 May 2022

Signature/Date