

Product Number:	102914
Lot Number:	210922
Expiration:	19-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Matt Smart

Name

QC

Title

MJA 3/15/2022

Signature/Date