

Product Number:	101476
Lot Number:	214903
Expiration:	29-May-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95 - 9.05

Review/Approval

Matt Smart

Name

QC

Title

MAT 06Jun2022

Signature/Date