

Product Number:	103847
Lot Number:	211313
Expiration:	23-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Matt Smart

Name

QC

Title

MRA 25M942022

Signature/Date