

Product Number:	103361
Lot Number:	215602
Expiration:	5-Jun-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Matt Smart

Name

QC

Title

Matt Smart 09/25/2022

Signature/Date