

Product Number:	101476
Lot Number:	308206
Expiration:	23-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95 - 9.05

Review/Approval

Matt Koulouris

Name

QC Supervisor

Title

[Signature] 29 Mar 2023

Signature/Date