

Product Number:	101556
Lot Number:	300607
Expiration:	6-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Matt Koulouris

QC Supervisor

[Signature] 17Jan2023

Name

Title

Signature/Date