

Product Number:	101474
Lot Number:	307613
Expiration:	17-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Matt Kodouris

Name

QC Supervisor

Title

MM 29 Mar 2023

Signature/Date