

Product Number:	104101
Lot Number:	220005
Expiration:	19-Jul-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Matt Koulouris
Name

QC Supervisor
Title

[Signature] 06Feb2023
Signature/Date