

Product Number:	101356
Lot Number:	227910
Expiration:	6-Oct-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

*Matt Koulouris*

Name

*QC Supervisor*

Title

*[Signature]* 10oct2022

Signature/Date