

Product Number:	101356
Lot Number:	227617
Expiration:	3-Oct-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Matt Koubaris

QC Supervisor

M. Koubaris 10 OCT 2022

Name

Title

Signature/Date