

Product Number:	101359
Lot Number:	227312
Expiration:	30-Sep-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Matt Koberis

Name

QC Supervisor

Title

[Signature] 10 Oct 2022

Signature/Date