

Product Number:	101359
Lot Number:	306906
Expiration:	10-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Matt Karburis

Name

QC Supervisor

Title

[Signature] 30 Mar 2023

Signature/Date