

Product Number:	101359
Lot Number:	223012
Expiration:	18-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Name

Title

Signature/Date

Sonia Khillan QC Manager Sonia Khillan 1/13SEP2022