

Product Number:	101353
Lot Number:	225116
Expiration:	8-Sep-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Matt Smart

Name

QC

Title

MAT 19sep2022

Signature/Date