

Product Number:	101359
Lot Number:	214208
Expiration:	22-May-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

Matt Smart 01/24/2022

Signature/Date