

Product Number:	101359
Lot Number:	217216
Expiration:	21-Jun-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

M. A. 27 Jun 2022

Signature/Date