

Product Number:	100453
Lot Number:	217503
Expiration:	24-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

Matt Smart
Name

QC

Title

MJA

27 Jun 2022

Signature/Date