

Product Number:	100453
Lot Number:	223714
Expiration:	25-Aug-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

Matt Smart

QC

MAT 19 Sep 2022

Name

Title

Signature/Date