

Product Number:	100376
Lot Number:	215209
Expiration:	1-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. Smart 06 Jun 2022

Signature/Date