

Product Number:	100458
Lot Number:	225608
Expiration:	13-Sep-2023
Storage:	2-8 °C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Matt Smart

Name

QC

Title

[Signature] 19Sep2022

Signature/Date