

Product Number:	101362
Lot Number:	211113
Expiration:	21-Apr-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

*Matt Smeit*

QC

*MJA 01 Jun 2022*

Name

Title

Signature/Date