

Product Number:	101954
Lot Number:	231416
Expiration:	10-Nov-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Killan
Name

QC Manager
Title

Sonia Killan / 14 NOV 2022
Signature/Date