

Product Number:	103158
Lot Number:	212413
Expiration:	4-May-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval Matt Smart QC MSA 10 May 2022
Name Title Signature/Date