

Product Number:	100376
Lot Number:	225011
Expiration:	7-Sep-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan QC Manager Sonia Khillan / 04 OCT 2022
Name Title Signature/Date