

Product Number:	101357
Lot Number:	303211
Expiration:	1-Feb-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 13 FEB 2023
Name Title Signature/Date