

Product Number:	100453
Lot Number:	233906
Expiration:	5-Dec-2023
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 12/21/2022
Name Title Signature/Date