

Product Number:	101352
Lot Number:	234312
Expiration:	9-Dec-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 103 JAN 2023
Name Title Signature/Date