

Product Number:	100458
Lot Number:	312905
Control Number:	312905A
Expiration:	9-May-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval Sonia Khillan QC Manager Sonia Khillan 23MAY2023
Name Title Signature/Date