

Product Number:	101475
Lot Number:	313205
Control Number:	313205A
Expiration:	12-May-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.40	7.35 - 7.45	PASS

Review/Approval Sonia Khillan GC Manager Soni K. Khillan 01SEP2023
 Name Title Signature/Date