





Product Number:	101475
Lot Number:	326106
Expiration:	18-Sep-2024
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Name: Prakash Jett EAC Prakash
 Title: SUPERVISOR
 Signature/Date: 27 SEP 2023