





Product Number:	101474
Lot Number:	326408
Expiration:	21-Sep-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.41	7.35 - 7.45	PASS

Review/Approval




 Name: Pratik Jetti Title: SUPERVISOR Signature/Date:  27 SEP 2023