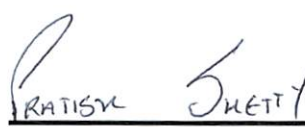




Product Number:	103847
Lot Number:	328318
Expiration:	10-Oct-2024
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval



 Name: Pratik Shetty Title: SUPERVISOR Signature/Date:  08 DEC 2023