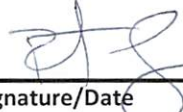


Product Number:	101357
Lot Number:	402306
Expiration:	23-Jul-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.19	6.18 - 6.22	PASS

Review/Approval

ROBERT SOMOGYI QC MANAGER  25 JAN 2024

---

Name Title Signature/Date