

Product Number:	100453
Lot Number:	513606
Expiration:	16-Nov-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Kathleen Corneli
Name
Quality Control
Supervisor
Title


18 Aug 2025
Signature/Date