

Product Number:	103359
Lot Number:	517101
Expiration:	20-Jun-2026
Storage:	2 - 8°C

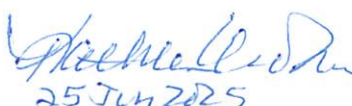
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.98	7.95 - 8.05	PASS

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


25 Jun 2025
Signature/Date