

Product Number:	103358
Lot Number:	517403
Expiration:	23-Jun-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.04	7.95 - 8.05	PASS

Review/Approval

Kathleen
Cornell

Name

Quality Control
Supervisor

Title


26 Jun 2025

Signature/Date