

Product Number:	103360
Lot Number:	517102
Expiration:	20-Jun-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.96	7.95 - 8.05	PASS

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

Kathleen Cornell
25 Jun 2025
Signature/Date