

Product Number:	102910
Lot Number:	523401
Expiration:	22-Aug-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.01	7.95 - 8.05	PASS

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

Kathleen Cornell
02 Sep 2025

Signature/Date