

Product Number:	102908
Lot Number:	523408
Expiration:	22-Aug-2026
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


04 Sep 2025
Signature/Date