

Product Number:	103158
Lot Number:	507621
Expiration:	17-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.35	7.35 - 7.45	PASS

Review/Approval

Heidi Caswell
Name

Director, QC
Title

Heidi Caswell 05 Jun 2025
Signature/Date