

Product Number:	101355
Lot Number:	509412
Expiration:	4-Oct-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature]
29 May 2025

Signature/Date