

Product Number:	101475
Lot Number:	521907
Expiration:	7-Feb-2027
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.40	7.35 - 7.45	PASS

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature] 18 Aug 2025

Signature/Date