

Product Number:	100457
Lot Number:	513908
Expiration:	19-Nov-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature] 29 May 2025

Signature/Date