

Product Number:	100457
Lot Number:	521020
Expiration:	29-Jan-2027
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]
19 Nov 2025

Signature/Date