

Product Number:	100458
Lot Number:	512007
Expiration:	30-Oct-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature] *23 May 2025*

Signature/Date