

Product Number:	100460
Lot Number:	510809
Expiration:	18-Oct-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature] 23 May 2025

Signature/Date