

Product Number:	100460
Lot Number:	510810
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

M. J. M. 18 Aug 2025
Signature/Date