

Product Number:	104101
Lot Number:	515315
Expiration:	2-Dec-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Heidi Caswell
Name

Director, QC
Title

Heidi Caswell 05 Jun 2025
Signature/Date