

Product Number:	101358
Lot Number:	523110
Expiration:	19-Feb-2027
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]
05 Nov 2025

Signature/Date