

Product Number:	101358
Lot Number:	507117
Expiration:	12-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi

QC Manager

 14 MAR 2025

Name

Title

Signature/Date