

Product Number:	101358
Lot Number:	507118
Expiration:	12-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

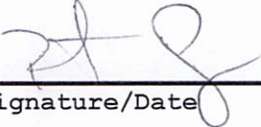
Robert Somogyi

QC Manager

Name

Title

Signature/Date

 07 APR 2025