

Product Number:	100458
Lot Number:	506316
Expiration:	4-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Robert Somogyi

QC Manager

 24 MAR 2025

Name

Title

Signature/Date