

Product Number:	100458
Lot Number:	508604
Expiration:	27-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

Name

Title

Signature/Date

 31 MAR 2025