

Product Number:	101955
Lot Number:	506314
Expiration:	4-Sep-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.43	7.35 - 7.45	PASS

Review/Approval

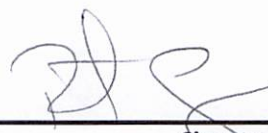
Robert Somogyi

QC Manager

Name

Title

Signature/Date

 18 MAR 2025