

Product Number:	103358
Lot Number:	506414
Expiration:	5-Mar-2026
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.96	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 10 MAR 2025

Name

Title

Signature/Date