

Product Number:	103158
Lot Number:	407912
Expiration:	19-Sep-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.38	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 25 MAR 2024

Signature/Date