

Product Number:	100457
Lot Number:	402510
Expiration:	25-Jul-2025
Storage:	2-8°C

**Laboratory Analysis**

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

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 02 FEB 2024

Signature/Date