

Product Number:	100457
Lot Number:	411728
Expiration:	26-Oct-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 01MAY2024

Signature/Date