

Product Number:	101472
Lot Number:	402613
Expiration:	26-Jul-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.39	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 09 FEB 2024

Signature/Date