

Product Number:	103847
Lot Number:	325907
Expiration:	16-Mar-2025
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 07 FEB 2024

Signature/Date