

Product Number:	103847
Lot Number:	432002
Expiration:	15-May-2026
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager

 21 NOV 2024

Name

Title

Signature/Date