

Product Number:	103847
Lot Number:	409214
Expiration:	1-Oct-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

QC Manager

Name

Title

 26 JUN 2024
Signature/Date