


Product Number:	101956
Lot Number:	404317
Expiration:	12-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.43	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI QC MANAGER
Name Title

 19 MAR 2024
Signature/Date