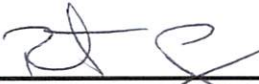


Product Number:	101955
Lot Number:	422015
Expiration:	7-Feb-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.43	7.30 - 7.50	PASS

Review/Approval

Robert Somogyi	QC Manager	 16 AUG 2024
Name	Title	Signature/Date