

Product Number:	102907
Lot Number:	504152
Expiration:	6-Feb-2025
Storage:	2-8°C

Component	Item Number	Lot Number
Lysate Diluent A	102908	404703
Lysate Diluent B	102910	403706
Lysate Diluent C	102912	403708
Lysate Diluent D	102914	403710

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title



Signature/Date

21 FEB 2024