

Product Number:	102907
Lot Number:	504479
Expiration:	2-Dec-2025
Storage:	2-8°C

Component	Item Number	Lot Number
Lysate Diluent A	102908	433704
Lysate Diluent B	102910	433719
Lysate Diluent C	102912	434008
Lysate Diluent D	102914	434010

Laboratory Analysis

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

Review/Approval

Kyle Giroux
Name

Senior Document Control Specialist, QA
Title

[Signature] 16 Dec 2024
Signature/Date