

Product Number:	102907
Lot Number:	504717
Expiration:	22-Aug-2026
Storage:	2-8°C

Component	Item Number	Lot Number
Lysate Diluent A	102908	523408
Lysate Diluent B	102910	523401
Lysate Diluent C	102912	523404
Lysate Diluent D	102914	523402

Laboratory Analysis

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

Review/Approval

Signature

DR / 09 Sep 2025

Name

DR

Title

QA Specialist II, QA