

Product Number:	102907
Lot Number:	504496
Expiration:	24-Jun-2025
Storage:	2-8°C

Component	Item Number	Lot Number
Lysate Diluent A	102908	417606
Lysate Diluent B	102910	417701
Lysate Diluent C	102912	417608
Lysate Diluent D	102914	434010

**Laboratory Analysis**

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

Review/Approval

Dhruvi Patel      Doc Control I, QA      SDP 23 Jan 2025

Name

Title

Signature/Date