

Product Number:	102907
Lot Number:	504267
Expiration:	3-May-2025
Storage:	2-8°C

Component	Item Number	Lot Number
Lysate Diluent A	102908	412401
Lysate Diluent B	102910	412403
Lysate Diluent C	102912	412405
Lysate Diluent D	102914	412407

Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager



16 MAY 2024

Name

Title

Signature/Date