

<b>Product Number:</b>	102907
<b>Lot Number:</b>	504305
<b>Expiration:</b>	3-May-2025
<b>Storage:</b>	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	412421

**Laboratory Analysis**

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

Review/Approval

Robert Somogyi

QC Manager

 02JUL2024

Name

Title

Signature/Date