

Product Number:	103360
Lot Number:	434009
Expiration:	5-Dec-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.04	7.95 - 8.05	PASS

Review/Approval

*Robert Somogyi*

QC MANAGER

*[Signature]* 16 Dec 2024

Name

Title

Signature/Date

# Certificate of Testing Review Checklist

Component Name:	Lysate Diluent C
Product Number:	103360
Lot Number:	434009

Content	Data Entered By (Initial/Date)	Data Checked By (Initial/Date)
		JV 12 DEC 2024
Product Name	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Product Number	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Lot Number	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Expiration	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Storage Condition	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Laboratory Analysis:		
Characteristic	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Means of Analysis	Accept <input type="checkbox"/> Reject <input type="checkbox"/> N/A <input checked="" type="checkbox"/>	Accept <input type="checkbox"/> Reject <input type="checkbox"/> N/A <input checked="" type="checkbox"/>
Specification	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Result	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>
Pass/Fail	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>	Accept <input checked="" type="checkbox"/> Reject <input type="checkbox"/> N/A <input type="checkbox"/>

MISSER Box PDS 16 DEC 2024

Comments: \_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_

N/A JV 12 DEC 2024

**1. REVISION HISTORY**

Revision	Detailed Description	Originator
1.0	Initial Release	D. Rice
02	DCR-16-0600: Transfer documents from QMS 2.0 to R&D vault. Archive the above docs in QMS 2.0 once released in R&D vault.	S. Chin
03	DCR-22-3411: Fixed macros in the header of the document. No change to content.	S. Gallagher
04	DCR-23-1983: - Updated heading title to "Certificate of Testing" - Removed "(Explain any N/A selections)"	V. Velasquez

*End of Document*