

Product Number:	102912
Lot Number:	207612
Expiration:	17-Mar-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia Khillan / 19 JUL 2022
Signature/Date

Component Certificate of Analysis Review Checklist

Component Name:	Lysate Diluent C (30 mL)
Product Number:	102912
Lot Number:	207612

Content	Data Entered By (Initial/Date)	Data Checked By (Initial/Date)
		WV / 14 JUL 2022
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"/>

Comments (Explain any N/A

selections):

~~N/A~~

~~19 JUL 2022 SK~~

REVISION HISTORY

Revision	Detailed Description	Date	Originator
1.0	Initial Release	17Dec2015	D. Rice
1.1	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.	28Sep2016	S. Chin

End of Document