

Beckman Coulter™ SYNCHRON CX® Systems

CAT NO:

APPLICATION PROCEDURE

TR66056

Reagent Preparation: Ready-to-use

Sample, Control, and Calibrator Preparation: An off-line dilution is required. All samples, controls, and calibrators should be prediluted 1:10 [sample to diluent ratio] with distilled or deionized water prior to assaying. For example, dilute 5 µL of sample with 50 µL of diluent.

INSTRUMENT PARAMETERS:

Test Name	LI
Reaction Type	ENDPOINT 2
Units	mmol/L
Precision	X.XX
Reaction Direction	Negative
Calculation Factor	0
Math Model	Linear
Calibration Time Limit	168 hr
Number of Calibrators	1
#1	1.00 [+]
Primary Wavelength	520 nm
Secondary Wavelength	470 nm
Sample Volume	3 µL

REAGENTS:

Primary Inject (first)/First Inject	
Compartment/Component	B
Volume/Dispense Volume	300 µL
Add Time/Inject Time	N/A
Primary Inject (second)/Second Inject	
Compartment/Component	None
Volume/Dispense Volume	0 µL
Add Time/Inject Time	N/A
Secondary Inject/Third Inject	
Compartment/Component	N/A
Volume/Dispense Volume	0 µL
Add Time/Inject Time	N/A
Blank	
Start Read	272 sec
End Read	304 sec
Reaction 1	
Start Read	128 sec
End Read	160 sec
Reaction 2	
Start Read	N/A
End Read	N/A

USABLE RESULT RANGE:

SYNCHRON CX	
Lower Limit	0.04
Upper Limit	3.00

ERROR DETECTION LIMITS:

Reagent Blank/Blank	
ABS Low Limit	0.000
ABS High Limit♦	1.500
Rate Low Limit	N/A
Rate High Limit	N/A
Mean Deviation	N/A
Reaction/Reaction 1	
ABS Low Limit	-0.700
ABS High Limit ♦	1.500
Rate Low Limit	N/A
Rate High Limit	N/A
Mean Deviation	N/A

SUBSTRATE DEPLETION:

Initial Rate♦	-99.999
Delta ABS♦	1.500

MULTI POINT SPAN: 1-2: -0.001

+ see recommended Lithium Standard TR66901 [1 mmol/L]

♦ Default parameter

***This is a preliminary instrument application and has not been validated or verified on the Beckman Coulter™ SYNCHRON CX® Systems analyzer. Each laboratory using this application is responsible for validating and verifying the performance of the application on their system.**

For further information on the use of this reagent, please consult the product insert, or contact Technical Support: U.S. 1-800-528-0494, Option 2, International 1-540-869-3200

Fisher Diagnostics, a division of Fisher Scientific, a part of Thermo Fisher Scientific, 8365 Valley Pike, Middletown, VA 22645 USA

© 2008 Thermo Fisher Scientific Inc. All rights reserved. SYNCHRON CX® is a trademark of Beckman Coulter Inc., Fullerton, CA 92835. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. www.thermo.com/diagnostics