

Beckman Coulter™ SYNCHRON CX® Systems

CAT NO:

APPLICATION PROCEDURE

TR25421

Reagent Preparation: Ready-to-use

INSTRUMENT PARAMETERS:

Test Name	AMY
Reaction Type	RATE 1
Units	U/L
Precision	X
Reaction Direction	POSITIVE
Calculation Factor	10933
Math Model	Linear
Calibration Time Limit	0
Number of Calibrators	0
Primary Wavelength	410 nm
Secondary Wavelength	520 nm
Sample Volume	6 µL

REAGENTS:

Primary Inject (first)/First Inject	
Compartment/Component	A
Volume/Dispense Volume	240 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	32 sec
End Read	64 sec
Reaction 1	
Start Read	64 sec
End Read	192 sec
Reaction 2	
Start Read	n/a
End Read	n/a

USABLE RESULT RANGE:

SYNCHRON CX	
Lower Limit	10
Upper Limit	2000

ERROR DETECTION LIMITS:

Reagent Blank/Blank	
ABS Low Limit	-0.010
ABS High Limit	0.300
Rate Low Limit	N/A
Rate High Limit	N/A
Mean Deviation	N/A
Reaction/Reaction 1	
ABS Low Limit	0.000
ABS High Limit♦	1.500
Rate Low Limit	N/A
Rate High Limit	N/A
Mean Deviation	N/A

SUBSTRATE DEPLETION:

Initial Rate	1.000
Delta ABS♦	1.500

MULTI POINT SPAN: n/a

♦ Default setting

***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