

**Beckman Coulter™ SYNCHRON LX®/Unicel® DxC 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	n/a
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	-80 sec
End Read	-20 sec
Reaction 1	
Start Read	60 sec
End Read	180 sec
Reaction 2	
Start Read	n/a
End Read	n/a

**USUABLE RESULT RANGE:**

Lower Limit	10
Upper Limit	2000

**ERROR DETECTION LIMITS:**

## Reagent Blank/Blank

ABS Low Limit ♦	-1.500
ABS High Limit♦	2.200
Rate Low Limit♦	-1.500
Rate High Limit♦	2.200
Mean Deviation♦	2.200

## Reaction/Reaction 1

ABS Low Limit ♦	-1.500
ABS High Limit ♦	2.200
Rate Low Limit ♦	-1.500
Rate High Limit ♦	2.200
Mean Deviation ♦	2.200

## Reaction/Reaction 2

ABS Low Limit ♦	n/a
ABS High Limit ♦	n/a
Rate Low Limit ♦	n/a
Rate High Limit ♦	n/a
Mean Deviation ♦	n/a

**SUBSTRATE DEPLETION:**

Initial Rate♦	99.999
Delta ABS♦	2.200

♦ Default settings

**\*This is a preliminary instrument application and has not been validated or verified on the Beckman Coulter™ SYNCHRON LX®/Unicel® DXC 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

© 2009 Thermo Fisher Scientific Inc. All rights reserved. SYNCHRON LX®/Unicel® 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](http://www.thermo.com/diagnostics)