

Olympus AU400® / AU600®

APPLICATION PROCEDURE

Reagent Preparation: Ready-to-use

CAT NO:
TR25421

Specific test parameters

Test No	#	Test name	AMY	Sample type	Serum	Page	1/2
Sample Volume	4	μL	Dil. Volume	0	μL	Min OD	Max OD
Reagent 1 Volume	160	μL	Dil. Volume	0	μL	L	H
Reagent 2 Volume		μL	Dil. Volume	0	μL	Reagent OD Limit	
Wavelength	Main	410	nm	Sub	520	nm	
Method					RATE		
Reaction					+		
Point 1	First	4		Last	10		
Point 2	First			Last			
Linearity	First	15	%	Second		%	
No Lag Time					YES		
						* Units = U/L	
						On Board Stability Period	#

Select using space key, or select from list displayed by guide key.

Specific test parameters

Test No	#	Test name	AMY	Sample type	Serum	Page	2/2
Value/Flag		Value		Level L	#	Level H	#
Normal range							
1 - 6	Sex	Age	L	Age	H	L	H
	#	# Y	# M	# Y	# M →	#	#
7	Non select					±	±
8	Out of range					#	#
Panic Value						L	H
						#	#

Select the function using the function key or the mouse.

Calibration Specific

Test No	#	Test name	AMY				
Cal type	1	MB		Count	2		
Formula	1	Y = AX + B		Process			
Selection calibrator							
Point 1	Cal No.	OD	Conc	Factor/OD-L	Factor/OD-H		
Point 2 - 7							
1 - Point cal. Point							
MB type factor		5313					
Calibrator Stability period		#					

Select the function using the function key of the mouse.

User defined parameters

± Normal Ranges - this parameter also sets decimal places for results

Olympus AU400® / AU600®
APPLICATION PROCEDURE

Reagent Preparation: Ready-to-use

CAT NO:
TR25421

***This is a preliminary instrument application and has not been validated or verified on the Olympus AU400®/AU600® 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. AU is a trademark of Olympus America, Inc. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. www.thermo.com/diagnostics