

Olympus AU400® / AU600®

APPLICATION PROCEDURE

Reagent Preparation: Ready-to-use

CAT NO:

TR66056

Specific test parameters										
Test No	#	Test name	LI	Sample type	Serum	Page	1/2			
Sample Volume	2	μL	Dil. Volume	0	μL	Min OD	Max OD			
Reagent 1 Volume	150	μL	Dil. Volume	0	μL	L		H		
Reagent 2 Volume	0	μL	Dil. Volume	0	μL	Reagent OD Limit				
Wavelength	Main	520	nm	Sub	480	nm	First L	-2.0000	First H	2.5000
		Last L	-2.0000		Last H	2.5000				
Method					END	Dynamic Range				
Reaction					-	L	0.04	H	3.00	
Point 1	First	0	Last	6	Correlation Factor				A	1.0
Point 2	First		Last						B	0.0
Linearity	First		%	Second		%	* Units = mmol/L		Sample Pre-dilution rate	10
No Lag Time						On Board Stability Period		#		

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

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

Select the function using the function key or the mouse.

Test No	#	Test name	LI						
Cal type	8	AB	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									
Calibrator Stability period		#							

Select the function using the function key of the mouse.

# User defined parameter

± Normal ranges - this parameter also set decimal places for results

Olympus AU400® / AU600®

**APPLICATION PROCEDURE**

**Reagent Preparation:** Ready-to-use

**CAT NO:**

TR66056

**\*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](http://www.thermo.com/diagnostics)