

CEDIA[®] CYCLOSPORINE PLUS (LOW RANGE) APPLICATION BECKMAN COULTER AU480[®]/AU680[®]/AU5800[®]

Beckman Coulter Reagent REF A31849

The CEDIA Cyclosporine PLUS assay is for the in vitro quantitative determination of cyclosporine in human whole blood using automated clinical chemistry analyzers as an aid in the management of cyclosporine therapy in kidney, liver, and heart transplants.

For In Vitro Diagnostic Use Only

Purpose

The information provided in this application sheet is intended as a supplement to the package insert. Refer to the package insert for information on intended use, reagent storage, reagent preparation, specimen collection, specimen preparation, specimen storage, quality control, and additional performance data.

Ordering Information

Item	Size	Beckman Coulter Reorder Number
CEDIA Cyclosporine PLUS Assay	R1 41 mL, R2 19 mL, Lysing Reagent 98 mL, Low Cal A 2.5 mL, Low Cal B 2.5 mL	A31849
More Diagnostics Rap/Tac/CsA Control Level 1	4 x 4 mL	B51007
More Diagnostics Rap/Tac/CsA Control Level 2	4 x 4 mL	A53712
More Diagnostics Rap/Tac/CsA Control Level 3	4 x 4 mL	A53713
AU Bottle	30 mL	63094
AU Bottle	60 mL	63093

Technical Support

For Technical Support, please contact your local Beckman Coulter Representative.

Reagent Storage

Refer to the package insert for information on reagent storage.

Continued on next page

Instructions For Use

Procedure for Analyzer

Refer to the operator's manuals for information on analyzer operation. Refer to the package insert for complete reagent preparation.

Prior to pouring into AU bottles, allow the reagent to equilibrate for 15 minutes at refrigerated temperature (2 to 8°C). Dispense R1 reagent and R2 reagent into appropriate AU bottles as shown in the table below:

CEDIA Cyclosporine Assay Kit	AU Reagent Bottle	
	R1 Compartment	R2 Compartment
Antibody/Substrate Reagent [R1]	One Bottle (60 mL)	
Enzyme Conjugate Reagent [R2]		One Bottle (30 mL)

Warning: These reagents have to be programmed to fixed positions. Do not use the Thermo reagent bottles directly on the AU analyzer.

If running both CsA Low Range & High Range, shared reagent can be set up as follows:

In the "Common Test Parameter" menu, select the "Test Name" tab. Enter the same Reagent ID for CSAL and CSAH in the Reagent ID column.

Results and Data Interpretation

Results for samples will be printed in ng/mL.

Specimen Preparation

Refer to the package insert for the complete specimen preparation. The product insert can be found at the Thermo Fisher website:

www.thermoscientific.com/Diagnostics

Calibration

Use the CEDIA Cyclosporine PLUS Low Range Calibrator Kit. The calibrators are prepared like patient samples. The value on the bottle is the value to use in the parameters below. These are lot number specific and should be updated when calibrator lot numbers change.

Application Parameters

Parameters The following tables outline the CEDIA Cyclosporine PLUS Assay chemistry parameters on the Beckman AU480, AU680, and AU5800 analyzers.

CEDIA CYCLOSPORINE PLUS – LOW RANGE, AU480

Specific Test Parameters																					
General		LIH		ISE		Range															
Test Name:		CSAL		<		>		Type:		Serum		Operation:		Yes							
Sample Volume		19.0		μL		Dilution		0		μL		OD Limit									
Pre-Dilution Rate		1								Min. OD		-2.00		Max. OD		3.00					
Reagents Volume: R1(R1-1)		146		μL		Dilution		0		μL		Reagent OD limit:									
										First Low		-2.00		High		3.00					
										Last Low		-2.00		High		3.00					
R2 (R2-1)		75		μL		Dilution		0		μL		Dynamic Range Low		25		High		450			
Wavelength: Pri.		570		nm		Sec.		660		nm		Correlation Factor A		1		B		0			
Method:		FIXED1		▽								Factor for Maker A		1		B		0			
Reaction slope:		+		▽								Onboard Stability		#		Days		#		Hour	
Measuring Point 1: First		24				Last		27				LIH Influence Check		#		▽					
Measuring Point 2: First						Last						Lipemia				▽					
Linearity:				%								Icterus				▽					
No Lag Time:		No		▽								Hemolysis				▽					

Specific Test Parameters															
General		ISE		Range											
Test Name:		CSAL		<		>		Type:		Serum					
Value/Flag:		#		▽		Level L:		#		Level H:		#			
Specific Ranges:															
		From		To		Low		High		Panic Value					
		Sex		Year		Month		Year		Month		Low		High	
<input type="checkbox"/> 1.		#		▽		#		#		#		#		#	
<input type="checkbox"/> 2.		#		▽		#		#		#		#		#	
<input type="checkbox"/> 3.		#		▽		#		#		#		#		#	
<input type="checkbox"/> 4.		#		▽		#		#		#		#		#	
<input type="checkbox"/> 5.		#		▽		#		#		#		#		#	
<input type="checkbox"/> 6.		#		▽		#		#		#		#		#	
<input type="checkbox"/> 7.		No demographics								#		#			
<input type="checkbox"/> 8.		Not within expected values								#		#			
Unit		ng/mL				Decimal Places		#							

Continued on next page

CEDIA CYCLOSPORINE PLUS – LOW RANGE, AU480, continued

Calibration Specific									
General		ISE							
Test Name:		CSAL	<	>	Type	Serum	<input type="checkbox"/> Use Serum Cal.		
Calibration Type:		AA	Formula:		Y = AX+B	Counts:	2		
<Calibrator Parameters>									
	Calibrator †	OD	Conc	Factor Range		Slope Check			
				Low	High				
Point 1:	#		*	-99999	99999	Allowable Range Check			
Point 2:	#		*			<input type="checkbox"/> Reagent Blank			
Point 3:						<input type="checkbox"/> Calibration			
Point 4:						Advanced Calibration			
Point 5:						Operation			
Point 6:						Interval (RB/ACAL)			
Point 7:									
Point 8:									
Point 9:									
Point 10:									
<Point Cal. For Master Curve>		No. of Correction Points			Use Master Curve	<input type="checkbox"/> Lot Calibration			
	Calibrator	OD	Conc	OD Range		Stability			
				Low	High	Reagent Blanks Calibration			
Point 1:						#	Day	#	Hour
Point 2:						#	Day	#	Hour
MB Type Factor:			1-Point Calibration Point			<input type="checkbox"/> With CONC-0			

User defined

* Lot specific calibrator values

Continued on next page

CEDIA CYCLOSPORINE PLUS – LOW RANGE, AU680

Specific Test Parameters																					
General		LIH		ISE		Range															
Test Name:		CSAL		<		>		Type:		Serum		Operation:		Yes							
Sample Volume		19.0		μL		Dilution		0		μL		OD Limit									
Pre-Dilution Rate		1								Min. OD		-2.00		Max. OD		3.00					
Reagents Volume:		R1(R1-1)		146		μL		Dilution		40		μL		Reagent OD limit:							
												First Low		-2.00		High		3.00			
														Last Low		-2.00		High		3.00	
R2 Volume		75		μL		Dilution		0		μL		Dynamic Range Low		25		High		450			
Common Reagent		Type		None		Name				Correlation Factor A		1		B		0					
Wavelength:		Pri.		570		nm		Sec.		660		nm		Factor for Maker A		1		B		0	
Method:		FIXED1																			
Reaction slope:		+								Onboard Stability		#		Days		#		Hour			
Measuring Point 1:		First		24		Last		27		LIH Influence Check		#									
Measuring Point 2:		First				Last				Lipemia											
Linearity:				%						Icterus											
No Lag Time:		No								Hemolysis											

Specific Test Parameters															
General		ISE		Range											
Test Name:		CSAL		<		>		Type:		Serum					
Value/Flag:		#		Level L:		#		Level H:		#					
Specific Ranges:															
		From		To		Low		High		Panic Value					
		Sex		Year		Month		Year		Month		Low		High	
<input type="checkbox"/> 1.		#		#		#		#		#		#		#	
<input type="checkbox"/> 2.		#		#		#		#		#		#		#	
<input type="checkbox"/> 3.		#		#		#		#		#		#		#	
<input type="checkbox"/> 4.		#		#		#		#		#		#		#	
<input type="checkbox"/> 5.		#		#		#		#		#		#		#	
<input type="checkbox"/> 6.		#		#		#		#		#		#		#	
		7. No demographics								#		#			
		8. Not within expected values								#		#			
Unit		ng/mL		Decimal Places		#									

Continued on next page

CEDIA CYCLOSPORINE PLUS – LOW RANGE, AU680, continued

Calibration Specific									
General		ISE							
Test Name:		CSAL		<	>	Type	Serum		<input type="checkbox"/> Use Serum Cal.
Calibration Type:		AA		Formula:		Y = AX+B		Counts:	2
<Calibrator Parameters>									
	Calibrator †	OD	Conc	Factor Range		Slope Check		+	
				Low	High			Allowable Range Check	
Point 1:	#		*	-99999	99999			<input type="checkbox"/> Reagent Blank	
Point 2:	#		*					<input type="checkbox"/> Calibration	
Point 3:								Advanced Calibration	
Point 4:								Operation	
Point 5:								No	
Point 6:								Interval (RB/ACAL)	
Point 7:									
Point 8:									
Point 9:									
Point 10:									
<Point Cal. For Master Curve>				No. of Correction Points		Use Master Curve		<input type="checkbox"/> Lot Calibration	
	Calibrator	OD	Conc	OD Range		Stability		Reagent Blanks Calibration	
				Low	High			# Day # Hour	
Point 1:								# Day # Hour	
Point 2:									
MB Type Factor:				1-Point Calibration Point				<input type="checkbox"/> With CONC-0	

User defined

* Lot specific calibrator values

Continued on next page

CEDIA CYCLOSPORINE PLUS – LOW RANGE, AU5800

Parameters		Specific Test Parameters									
General	LIH	ISE	HbA1c		Calculated Test	Range					
Test Name:		CSAL	<	>	Type:	Serum	Operation	Yes			
Sample Volume		17	μL	Dilution	0	μL	OD Limit				
Pre-Dilution Rate		1	∇	Diluent Bottle	#	∇	Min.OD	-2.00	Max.OD	3.00	
Rgt. Volume	R1(R1-1)	131	μL	Dilution	0	μL	Reagent OD Limit				
	R1-2		μL	Dilution		μL	First	Low	-2.00	High	3.00
							Last	Low	-2.00	High	3.00
	R2(R2-1)	67	μL	Dilution	0	μL	Dynamic Range Low		25	High	450
Common Rgt. Type		None		Name			Correlation Factor A	1	B	0	
Wavelength	Pri	570	∇nm	Sec.	660	∇nm	Factor for Maker A	1	B	0	
Method		FIXED1	∇				Onboard Stability Period	#	Day	#	Hour
Reaction Slope		+	∇				LIH Influence Check	#	∇		
Measuring Point1 1 st		24		Last	27		Lipemia		∇		
Measuring Point2 1 st				Last			Icterus		∇		
Linearity Limit			%				Hemolysis		∇		
Lag Time Check		No	∇								

Parameters		Specific Test Parameters						
General	LIH	ISE	HbA1c		Calculated Test	Range		
Test Name:		CSAL	<	>	Type:	Serum		
Value/Flag:		#	∇					
Specific Ranges:		From		Level To		Low	High	
	Sex	Year	Month	Year	Month	Low	High	
o 1.	#	∇	#	#	#	#	#	
o 2.	#	∇	#	#	#	#	#	
o 3.	#	∇	#	#	#	#	#	
o 4.	#	∇	#	#	#	#	#	
o 5.	#	∇	#	#	#	#	#	
o 6.	#	∇	#	#	#	#	#	
7.	Standard demographics					#	#	
8.	Not within expected values					#	#	
Panic Value	Low	#	High	#	Unit	ng/mL	Decimal Places	#

Continued on next page

CEDIA CYCLOSPORINE PLUS – LOW RANGE, AU5800, continued

Parameters		Calibration Parameters					
Calibrators		Calibration Specific					
General		ISE					
Test Name:		CSAL	<	>	Type: Serum	Cuvette .	
		<input type="checkbox"/> Use Serum Cal.					
Calibration Type:		AA	Formula: Y=AX+B		Counts: 2		
<Calibrator Parameters>		Range					
	Calibrator	OD	Conc	Low	High	Slope Check	
Point 1:	#		*	-99999	99999	<input type="text" value="+"/>	
Point 2:	#		*			Allowance Range Check	
Point 3:						<input type="checkbox"/> Reagent Blank	
Point 4:						<input type="checkbox"/> Calibration	
Point 5:						Advanced Calibration	
Point 6:						Operation	
Point 7:						<input type="text" value="No"/>	
Point 8:						Interval (RB/ACAL)	
Point 9:						<input type="text"/>	
Point 10:						<input type="text"/>	
<Point Cal. For		No. of Correction Points		Use Master Curve		<input type="checkbox"/> Lot Calibration	
Master Curve>		OD Range					
	Calibrator	OD	Conc	Low	High	Stability	
Point-1						Reagent Blank	# Day # Hour
Point-2						Calibration	# Day # Hour
MB Type Factor:		1-Point Calibration Point		None		<input type="checkbox"/> with Conc-0	

User defined

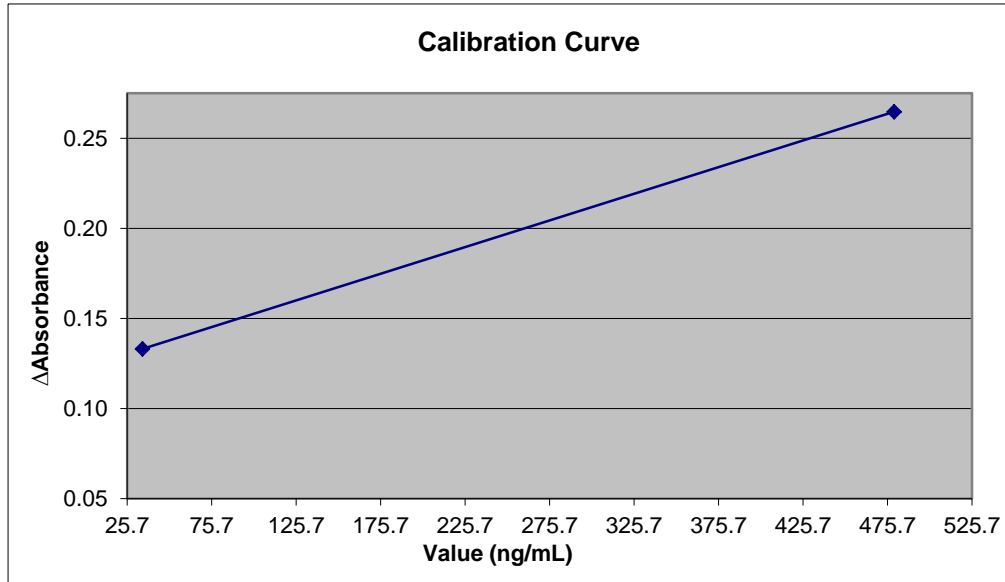
* Lot specific calibrator values

Results and Data Interpretation

**Performance
Data**

Refer to the CEDIA Cyclosporine assay kit package insert for additional information on results and data interpretation.

Example Calibration Curve, CsA LR (AU480):



Continued on next page

Precision

These degrees of precision and equivalency were obtained in typical testing procedures on an AU system and are not intended to represent the performance specifications for this reagent.

Control samples were tested in replicates of 2, twice per day for 20 days, total N = 80. The results are presented in the following table:

Controls	Control 1	Control 2	Control 3
AU480			
Mean (ng/mL)	77.2	220.6	347.5
Within-Run SD (ng/mL)	2.41	3.02	3.27
Within-Run CV (%)	3.1	1.4	0.9
Total SD (ng/mL)	6.03	11.45	16.37
Total CV (%)	7.8	5.2	4.7
AU680			
Mean (ng/mL)	48.0	201.7	309.4
Within-Run SD (ng/mL)	3.40	4.54	5.36
Within-Run CV (%)	7.1	2.3	1.7
Total SD (ng/mL)	6.12	7.69	10.36
Total CV (%)	12.7	3.8	3.3
AU5800			
Mean (ng/mL)	79.1	222.8	349.7
Within-Run SD (ng/mL)	3.51	7.03	5.61
Within-Run CV (%)	4.4	3.2	1.6
Total SD (ng/mL)	6.10	12.79	18.57
Total CV (%)	7.7	5.7	5.3

Continued on next page

Linearity

Ten levels of manufacturing calibrators were run against a single calibration curve and the linearity calculated for the AU480, AU680, and AU5800. The analytical range for this assay is 25 – 450 ng/mL. Error flags will appear for samples recovering above or below the assay range.

The Cyclosporine PLUS Low Range assay recovered between 102 – 105% of expected values on the AU480.

The Cyclosporine PLUS Low Range assay recovered between 95 – 100% of expected values on the AU680.

The Cyclosporine PLUS Low Range assay recovered between 91 – 98% of expected values on the AU5800.

LDD

A negative blood sample was run against the same calibration curve for 21 replicates. The LDD is calculated as $2 \times SD$.

The observed LDD for the Cyclosporine PLUS Low Range Assay was 0.01 ng/mL on the AU480.

The observed LDD for the Cyclosporine PLUS Low Range Assay was 0.04 ng/mL on the AU680.

The observed LDD for the Cyclosporine PLUS Low Range Assay was 0.01 ng/mL on the AU5800.

Accuracy and Correlation

One hundred and twelve blood samples were assayed with the CEDIA Cyclosporine PLUS Low Range Assay on the Beckman Coulter AU480 and tested with reference method Hitachi 911.

One hundred blood samples were assayed with the CEDIA Cyclosporine PLUS Low Range Assay on the Beckman Coulter AU680 and tested with reference method Hitachi 911.

One hundred and twelve blood samples were assayed with the CEDIA Cyclosporine PLUS Low Range Assay on the Beckman Coulter AU5800 and tested with reference method Hitachi 911.

A Deming's Regression Analysis for CsA LR yielded the following:

Beckman Coulter AU480 = $1.03 \times (\text{Hitachi 911}) - 0.40$ with a correlation coefficient of 0.995.

Beckman Coulter AU680 = $0.97 \times (\text{Hitachi 911}) + 13.00$ with a correlation coefficient of 0.995.

Beckman Coulter AU5800 = $1.00 \times (\text{Hitachi 911}) - 0.50$ with a correlation coefficient of 0.994.

Continued on next page

Additional Information

Important


Since Beckman Coulter does not manufacture the reagent or perform quality control or other tests on individual lots, Beckman Coulter cannot be responsible for the quality of the data obtained which is caused by performance of the reagent, any variation between lots of reagent, or protocol changes by the Manufacturer.

**Shipping
Damage**

Please notify your Beckman Coulter Clinical Support Center if this product is received damaged.

© 2016 Thermo Fisher Scientific, Inc. All rights reserved.
AU Series Systems are the registered trademarks of Beckman Coulter.
CEDIA® is a registered trademark of Roche.
All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries.



 Thermo Fisher Scientific Oy, Ratastie 2, P.O. Box 100, 01621 Vantaa, Finland
Tel: +358-9-329100/Fax: +358-9-32910300

End