

CEDIA[®] MYCOPHENOLIC ACID APPLICATION

Beckman Coulter DxC 700 AU[®]



Beckman Coulter Reagent REF B01460

Intended for the quantitative measurement of mycophenolic acid in human plasma using automated clinical chemistry analyzers as an aid in the management of mycophenolic acid therapy in renal and cardiac transplant patients.

For In Vitro Diagnostic Use Only
Rx 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	Catalog Number
CEDIA Mycophenolic Acid Assay	R1: 1 x 26 mL R2: 1 x 11 mL	B01460
CEDIA Mycophenolic Acid Calibrator Set	2 levels, 2 x 5 mL	B37609
MAS [®] Mycophenolic Acid Control 1 Kit	1 level, 4 x 5 mL	B37611
MAS Mycophenolic Acid Control 2 Kit	1 level, 4 x 5 mL	B01543
MAS Mycophenolic Acid Control 3 Kit	1 level, 4 x 5 mL	B01544
AU Bottle	20 X 15 mL	63165
AU Bottle	20 X 30 mL	63094

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 Mycophenolic Acid Assay Kit	AU Reagent Bottle	
	R1 Compartment	R2 Compartment
Antibody/Substrate Reagent [R1]	One Bottle (30 mL)	
Enzyme Conjugate Reagent [R2]		One Bottle (15 mL)

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

Results and Data Interpretation

Results for samples will be printed in µg/mL.

Specimen Preparation

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

www.thermofisher.com

Calibration

Use the CEDIA Mycophenolic Acid Calibrator set. The calibrators are prepared like patient samples. Refer to the value assignment card for calibrator values to program in the parameters below. These are lot number-specific and should be updated when calibrator lot numbers change.

CEDIA Mycophenolic Acid Assay Beckman Coulter System Parameters, DxC 700 AU

General		LIH		ISE		Calculated Test		Range	
Test Name:		#		Type:		Serum		Operation	
		▼				▼		Yes ▼	
Sample Volume	<input type="text" value="7.5"/>	μL	Dilution	<input type="text" value="0"/>	μL	OD Limit			
Pre-Dilution Rate	<input type="text" value="1"/>	▼			Min. OD	<input type="text" value="-2.0000"/>	Max OD	<input type="text" value="3.0000"/>	
Reagent Volume	R1 (R1-1)	<input type="text" value="150"/>	μL	Dilution	<input type="text" value="0"/>	μL	Reagent OD Limit	1 st	
	R1-2	<input type="text"/>	μL	Dilution	<input type="text"/>	μL	Last	Low	<input type="text" value="-2.0000"/>
	R2 (R2-1)	<input type="text" value="60"/>	μL	Dilution	<input type="text" value="0"/>	μL	Analytical Measuring Range	Low	<input type="text" value="0.30"/>
Common Reagent	Type	<input type="text" value="None"/>		Name	<input type="text" value="None"/>		Correlation Factor	A	<input type="text" value="1"/>
							B	<input type="text" value="0"/>	
Wavelength	Pri	<input type="text" value="570"/>	nm	Sec	<input type="text" value="660"/>	nm	Manufacturer Factor	A	<input type="text" value="1"/>
							B	<input type="text" value="0"/>	
Method	<input type="text" value="FIXED1"/>								
Reaction Slope	<input type="text" value="+"/>		▼		Onboard Stability Period	<input type="text" value="30"/>	Day	<input type="text" value="0"/>	Hour
Measuring Point-1	1st	<input type="text" value="24"/>	Last	<input type="text" value="27"/>	LIH Influence Check	<input type="text" value="No"/>			
Measuring Point-2	1st	<input type="text"/>	Last	<input type="text"/>	Lipemia	<input type="text" value="+"/>			
Linearity Limit	<input type="text"/>		%		Icterus	<input type="text" value="+"/>			
Lag Time Check	<input type="text" value="No"/>		▼		Hemolysis	<input type="text" value="+"/>			

General		LIH		ISE		Calculated Test		Range		
Test Name:		#		Type:		Serum		▼		
		▼				▼				
Value/Flag	<input type="text" value="#"/>		Level	Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>			
Specific Ranges										
	Sex	Year	From	Month	To	Year	Month	Other Type	Low	High
<input type="checkbox"/> 1:	<input type="text" value="#"/>	▼	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
<input type="checkbox"/> 2:	<input type="text" value="#"/>	▼	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
<input type="checkbox"/> 3:	<input type="text" value="#"/>	▼	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
<input type="checkbox"/> 4:	<input type="text" value="#"/>	▼	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
<input type="checkbox"/> 5:	<input type="text" value="#"/>	▼	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
<input type="checkbox"/> 6:	<input type="text" value="#"/>	▼	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
7:	Standard demographics								<input type="text" value="#"/>	<input type="text" value="#"/>
8:	Not within expected values								<input type="text" value="#"/>	<input type="text" value="#"/>
Critical Limits	Low	<input type="text" value="#"/>		High	<input type="text" value="#"/>		Unit	<input type="text" value="μg/mL"/>		
							Select	Decimal Places	<input type="text" value="2"/>	

User defined

CEDIA Mycophenolic Acid Assay Beckman Coulter System Parameters, DxC 700 AU, *continued*

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

User defined
* Lot specific calibrator values

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 Technical Support Center if this product is received damaged.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AU Series Systems are trademarks of Beckman Coulter.



EC REP B-R-A-H-M-S GmbH, Neuendorfstrasse 25, 16761, Hennigsdorf, Germany

End