

QMS[®] EVEROLIMUS APPLICATION Beckman Coulter AU680[®]/AU5800[®]

Beckman Coulter Reagent REF A53729 (US)

The QMS Everolimus assay is intended for the quantitative determination of everolimus in human whole blood on automated clinical chemistry analyzers. The results obtained are used as an aid in the management of kidney and liver transplant patients receiving everolimus therapy. This is an in vitro diagnostic device intended for clinical laboratory use only.

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 (US)
QMS [®] Everolimus Assay	R1 22 mL, R2 8 mL, Precipitation Reagent 8 mL	A53729
QMS Everolimus Calibrator Set	6 levels, 3mL 1 bottle ea	A53721
QMS Everolimus Control Set	3 levels, 3mL 1 bottle ea	A53717
AU Bottle	15 mL	63165
AU Bottle	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. Before use, invert several times, avoiding the formation of bubbles. Dispense R1 reagent and R2 reagent into appropriate AU bottles as shown in the table below:

	AU Reagent Bottle	
	R1 Compartment	R2 Compartment
QMS Everolimus Assay Kit		
Anti-Everolimus Polyclonal Antibody R1	One Bottle (30 mL)	
Everolimus-Coated Microparticles R2		One Bottle (15 mL)

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

Significant interference from QMS Everolimus into the Microalbumin (OSR6167) and Urine/CSF Albumin (B38858/B46435) assays has been observed due to reagent carryover in random access analyzers. Setup the recommended contamination parameters below:

AU680 Contamination Parameters							
No.	PRECEDING TEST NAME	FOLLOWING TEST NAME	REAGENT PROBE CLEANER KIND	WASH COUNT	EFFECTIVE OF WATER CLEANING	SAME USE	
						MIXER	CUVETTE
1	EVR	MALB/UALB	Water	3	Yes	Yes	No

AU5800 Contamination Parameters							
No.	PRECEDING TEST NAME	FOLLOWING TEST NAME	REAGENT PROBE CLEANER KIND	WASH COUNT	EFFECTIVE OF WATER CLEANING	SAME USE	
						MIXER	CUVETTE
1	EVR	MALB/UALB	Water	3	Yes	Yes	No

Note: For the AU5800 it is recommended to separate MALB and UALB from EVR by using designated rings, if possible.

**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. Due to sample stability, it is recommended not to exceed a maximum of 24 extracted samples per run. The product insert can be found at the Thermo Fisher website:

www.thermoscientific.com/Diagnostics

Calibration

Use the QMS Everolimus Calibrator Set. The calibrators are prepared like patient samples. The value on the bottle is the value to use in the parameters below. Lot-to-lot values do not change.

Application Parameters

Parameters

The following tables outline the QMS Everolimus Assay chemistry parameters on the Beckman AU680, and AU5800 analyzers.

QMS EVEROLIMUS, AU680

Specific Test Parameters									
General		LIH		ISE		Range			
Test Name:		EVER ▾		< >		Type: Serum ▾		Operation: Yes ▾	
Sample Volume		10.0 μL		Dilution 0 μL		OD Limit			
Pre-Dilution Rate		1				Min. OD -2.00		Max. OD 3.00	
Reagents Volume: R1(R1-1)		175 μL		Dilution 0 μL		Reagent OD limit:			
						First Low -2.00		High 3.00	
						Last Low -2.00		High 3.00	
R2 Volume		45 μL		Dilution 0 μL		Dynamic Range Low 2.0		High 20.0	
Common Reagent		Type None		Name		Correlation Factor A 1		B 0	
Wavelength:		Pri. 700 nm		Sec. None 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:		EVER ▾		< >		Type: Serum ▾			
Value/Flag:		# ▾		Level L: #		Level H: #			
Specific Ranges:									
		From		To		Low		High	
<input type="checkbox"/>	1.	Sex # ▾	Year #	Month #	Year #	Month #	#	#	#
<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		#			
Panic Value									
		Low #		High #					

Continued on next page

QMS EVEROLIMUS, AU680, continued

Calibration Specific									
General		ISE							
Test Name:		EVER	<	>	Type	Serum	<input type="checkbox"/> Use Serum Cal.		
Calibration Type:		6AB	Formula:		EIA Type 1	Counts:		#	
<Calibrator Parameters>									
	Calibrator †	OD	Conc	Factor Range		Slope Check			
				Low	High				
Point 1:	#		0.00	-2.0	3.0	Allowable Range Check			
Point 2:	#		1.50	-2.0	3.0	<input type="checkbox"/> Reagent Blank			
Point 3:	#		3.00	-2.0	3.0	<input type="checkbox"/> Calibration			
Point 4:	#		6.00	-2.0	3.0	Advanced Calibration			
Point 5:	#		12.00	-2.0	3.0	Operation			
Point 6:	#		20.00	-2.0	3.0	Interval (RB/ACAL)			
Point 7:									
Point 8:									
Point 9:									
Point10:									
<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			
Point 1:						Day			
Point 2:						Day			
MB Type Factor:		1-Point Calibration Point		<input type="checkbox"/> With CONC-0					

User defined

Continued on next page

QMS EVEROLIMUS, AU5800

Parameters		Specific Test Parameters									
General	LIH	ISE	HbA1c		Calculated Test	Range					
Test Name:		EVER	<	>	Type:	Serum	Operation	Yes			
Sample Volume	8.9	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)	156	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)	40	μL	Dilution	0	μL					
Common Rgt. Type	Pri	None	Name			Dynamic Range Low	2.0	High	20.0		
Wavelength		700	nm	Sec.	None	nm	Correlation Factor A	1	B	0	
Method		FIXED1					Factor for Maker A	1	B	0	
Reaction Slope		+				Onboard Stability Period	#	Day	#	Hour	
Measuring Point1 1 st		24		Last	27	LIH Influence Check	#				
Measuring Point2 1 st				Last		Lipemia					
Linearity Limit			%			Icterus					
Lag Time Check		No				Hemolysis					

Parameters		Specific Test Parameters						
General	LIH	ISE	HbA1c		Calculated Test	Range		
Test Name:		EVER	<	>	Type:	Serum		
Value/Flag:	#							
Specific Ranges:			Level		Low	#	High	#
	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.	Standard demographics						#	#
8.	Not within expected values						#	#
Panic Value	Low	#	High	#	Unit	ng/mL	Decimal Places	#

Continued on next page

QMS EVEROLIMUS, AU5800, continued

Parameters		Calibration Parameters							
Calibrators		Calibration Specific							
General		ISE							
Test Name:		EVER	<	>	Type	Serum	Cuvette .		
		<input type="checkbox"/> Use Serum Cal.							
Calibration Type:		6AB	Formula:		EIA Type 1	Counts:		2	
<Calibrator Parameters>		Range							
	Calibrator	OD	Conc	Low	High	Slope Check			
Point 1:	#		0.00	-2.0	3.0	-			
Point 2:	#		1.50	-2.0	3.0	Allowance Range Check			
Point 3:	#		3.00	-2.0	3.0	<input type="checkbox"/> Reagent Blank			
Point 4:	#		6.00	-2.0	3.0	<input type="checkbox"/> Calibration			
Point 5:	#		12.00	-2.0	3.0	Advanced Calibration			
Point 6:	#		20.00	-2.0	3.0	Operation			
Point 7:						No			
Point 8:						Interval (RB/ACAL)			
Point 9:									
Point 10:									
<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			
Point-2						Calibration			
MB Type Factor:		1-Point Calibration Point		None		<input type="checkbox"/> with Conc-0			

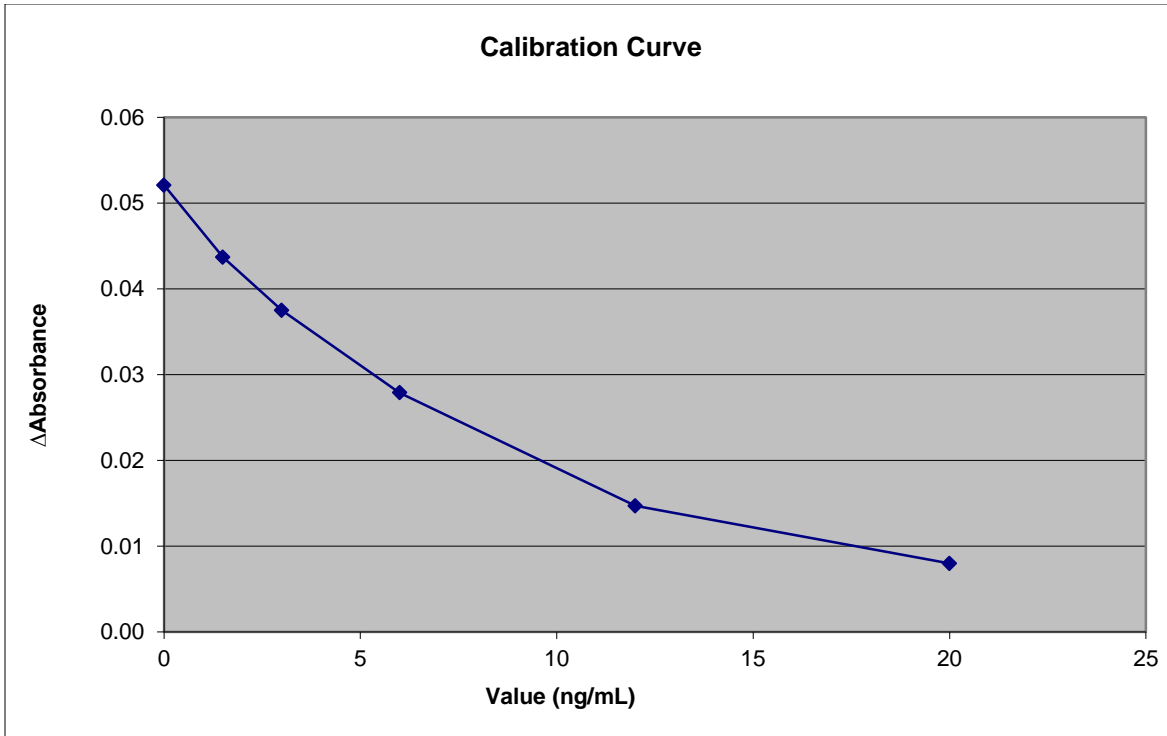
User defined

Results and Data Interpretation

Performance Data

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

Example Calibration Curve, Everolimus (AU680):



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
AU680			
Mean (ng/mL)	4.2	8.0	15.6
Within-Run SD (ng/mL)	0.17	0.35	0.81
Within-Run CV (%)	4.1	4.4	5.2
Total SD (ng/mL)	0.20	0.43	0.91
Total CV (%)	4.9	5.3	5.8
AU5800			
Mean (ng/mL)	3.9	7.4	14.1
Within-Run SD (ng/mL)	0.27	0.41	0.90
Within-Run CV (%)	6.8	5.5	6.4
Total SD (ng/mL)	0.40	0.64	1.14
Total CV (%)	10.1	8.7	8.1

Linearity

Eleven levels of calibrators and calibrator blends were run against a single calibration curve and the linearity calculated. The analytical range for this assay is 2.0 to 20.0 ng/mL. Error flags will appear for samples recovering above or below the assay range.

The Everolimus assay recovered between 97 – 106% of expected values on the AU680.

The Everolimus assay recovered between 93 – 103% of expected values on the AU5800.

Continued on next page

LDD

The negative calibrator was run against the same calibration curve for 21 replicates. The LDD is calculated as $2 \times \text{SD}$.

The observed LDD for the Everolimus Assay on the AU680 was 0.2 ng/mL.

The observed LDD for the Everolimus Assay on the AU5800 was 0.2 ng/mL.

Accuracy and Correlation

One hundred and fifty blood samples were assayed with the QMS Everolimus Assay on the Beckman Coulter AU680 and tested with reference method Hitachi 917.

One hundred and six blood samples were assayed with the QMS Everolimus Assay on the Beckman Coulter AU5800 and tested with reference method Hitachi 917.

A Deming's Regression Analysis for Everolimus yielded the following:

Beckman Coulter AU680 = $0.99 \times (\text{Hitachi 917}) - 0.13$ with a correlation coefficient of 0.923

Beckman Coulter AU5800 = $0.99 \times (\text{Hitachi 917}) + 0.13$ with a correlation coefficient of 0.982

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.

© 2014 Thermo Fisher Scientific, Inc. All rights reserved.
AU Series Systems are the registered trademarks of Beckman Coulter.
All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries.

End