

## QMS<sup>®</sup> EVEROLIMUS APPLICATION BECKMAN COULTER AU480<sup>®</sup> /AU680<sup>®</sup> /AU5800<sup>®</sup>

Beckman Coulter Reagent REF A53716 (International)

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 organ transplant patients receiving everolimus therapy.

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
QMS <sup>®</sup> Everolimus Assay	R1 22 mL, R2 8 mL, Precipitation Reagent 8 mL	A53716
QMS Everolimus Calibrator Set	6 levels, 3mL 1 bottle ea	A53724
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:

QMS Everolimus Assay Kit	AU Reagent Bottle	
	R1 Compartment	R2 Compartment
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 the 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:

AU480 Contamination Parameters									
No.	PRECEDING TEST NAME	Type	FOLLOWING TEST NAME	Type	REAGENT PROBE CLEANER KIND	WASH COUNT	EFFECTIVE OF WATER CLEANING	SAME USE	
								MIXER	CUVETTE
1	EVR	R1	MALB/UALB	R1	Water	3	Yes	Yes	No
2	EVR	R1	MALB/UALB	R2	Water	3	Yes	Yes	No

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](http://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 AU480, AU680, and AU5800 analyzers.

### QMS EVEROLIMUS, AU480

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 (R2-1)	45	μL	Dilution	0	μL	Dynamic Range Low	1.5	High	20.0		
Wavelength:	Pri.	700	nm	Sec.	None	nm	Correlation Factor A	1	B	0	
Method:	FIXED1										
Reaction slope:	+										
Measuring Point 1:	First	24	Last	27		Onboard Stability	#	Days	#	Hour	
Measuring Point 2:	First		Last			LIH Influence Check	#				
Linearity:											
No Lag Time:	No										
						Lipemia					
						Icterus					
						Hemolysis					

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

Continued on next page



**QMS EVEROLIMUS, AU480, continued**

Calibration Specific									
General		ISE							
Test Name:		EVER		<	>	Type	Serum		<input type="checkbox"/> Use Serum Cal.
Calibration Type:		6AB		Formula:		EIA Type 1		Counts:	2
<Calibrator Parameters>									
	Calibrator ↑	OD	Conc	Factor Range		Slope Check		-	
				Low	High			Allowable Range Check	
Point 1:	#		0.00	-2.0	3.0			<input type="checkbox"/> Reagent Blank	
Point 2:	#		1.50	-2.0	3.0			<input type="checkbox"/> Calibration	
Point 3:	#		3.00	-2.0	3.0			Advanced Calibration	
Point 4:	#		6.00	-2.0	3.0			Operation	
Point 5:	#		12.00	-2.0	3.0			Interval (RB/ACAL)	
Point 6:	#		20.00	-2.0	3.0				
Point 7:									
Point 8:									
Point 9:									
Point 10:									
<Point Cal. For Master Curve>									
	Calibrator	OD	Conc	OD Range		Use Master Curve		<input type="checkbox"/> Lot Calibration	
				Low	High			Stability	
Point 1:								Reagent Blanks	
Point 2:								Calibration	
								#	Day
								#	Hour
								#	Day
								#	Hour
MB Type Factor:		1-Point Calibration Point				<input type="checkbox"/> With CONC-0			

# User defined

*Continued on next page*



**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 1.5 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:	<table border="1"> <thead> <tr> <th></th> <th>From</th> <th>To</th> <th>Low</th> <th>High</th> </tr> <tr> <th></th> <th>Sex</th> <th>Year</th> <th>Month</th> <th>Year</th> <th>Month</th> <th>Low</th> <th>High</th> </tr> </thead> <tbody> <tr> <td><input type="checkbox"/> 1.</td> <td># ▾</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> </tr> <tr> <td><input type="checkbox"/> 2.</td> <td># ▾</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> </tr> <tr> <td><input type="checkbox"/> 3.</td> <td># ▾</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> </tr> <tr> <td><input type="checkbox"/> 4.</td> <td># ▾</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> </tr> <tr> <td><input type="checkbox"/> 5.</td> <td># ▾</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> </tr> <tr> <td><input type="checkbox"/> 6.</td> <td># ▾</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> <td>#</td> </tr> <tr> <td>7. No demographics</td> <td></td> <td></td> <td></td> <td></td> <td></td> <td>#</td> <td>#</td> </tr> <tr> <td>8. Not within expected values</td> <td></td> <td></td> <td></td> <td></td> <td></td> <td>#</td> <td>#</td> </tr> </tbody> </table>			From	To	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. No demographics						#	#	8. Not within expected values						#	#	<table border="1"> <thead> <tr> <th colspan="2">Panic Value</th> </tr> <tr> <th>Low</th> <th>High</th> </tr> </thead> <tbody> <tr> <td>#</td> <td>#</td> </tr> </tbody> </table>	Panic Value		Low	High	#	#
	From	To	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. No demographics						#	#																																																																															
8. Not within expected values						#	#																																																																															
Panic Value																																																																																						
Low	High																																																																																					
#	#																																																																																					
Unit	ng/mL	Decimal Places	#																																																																																			

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			Allowable Range Check	
Point 1:	#		0.00	-2.0	3.0			<input type="checkbox"/> Reagent Blank	
Point 2:	#		1.50	-2.0	3.0			<input type="checkbox"/> Calibration	
Point 3:	#		3.00	-2.0	3.0			Advanced Calibration	
Point 4:	#		6.00	-2.0	3.0			Operation	
Point 5:	#		12.00	-2.0	3.0			Interval (RB/ACAL)	
Point 6:	#		20.00	-2.0	3.0				
Point 7:									
Point 8:									
Point 9:									
Point 10:									
<Point Cal. For Master Curve>									
	Calibrator	OD	Conc	OD Range		Use Master Curve		<input type="checkbox"/> Lot Calibration	
				Low	High			Stability	
Point 1:								Reagent Blanks	
Point 2:								Calibration	
	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	None		Name			Dynamic Range Low	1.5	High	20.0		
Wavelength	Pri	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 <sup>st</sup>	24		Last	27				LIH Influence Check	#	▽	
Measuring Point2 1 <sup>st</sup>			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:	From	Level To		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: <input type="text" value="EVER"/> < > Type <input type="text" value="Serum"/> Cuvette . <input type="text"/>																																																																																																																													
<input type="checkbox"/> Use Serum Cal.																																																																																																																													
<table border="1" style="width:100%; border-collapse: collapse;"> <tr> <td colspan="2">Calibration Type: <input type="text" value="6AB"/></td> <td colspan="2">Formula: <input type="text" value="EIA Type 1"/></td> <td colspan="2">Counts: <input type="text" value="2"/></td> </tr> <tr> <td colspan="6" style="text-align: center;">&lt;Calibrator Parameters&gt;</td> </tr> <tr> <td></td> <td>Calibrator</td> <td>OD</td> <td>Conc</td> <td colspan="2" style="text-align: center;">Range</td> </tr> <tr> <td></td> <td></td> <td></td> <td></td> <td>Low</td> <td>High</td> </tr> <tr> <td>Point 1:</td> <td><input type="text" value="#"/></td> <td><input type="text"/></td> <td><input type="text" value="0.00"/></td> <td><input type="text" value="-2.0"/></td> <td><input type="text" value="3.0"/></td> </tr> <tr> <td>Point 2:</td> <td><input type="text" value="#"/></td> <td><input type="text"/></td> <td><input type="text" value="1.50"/></td> <td><input type="text" value="-2.0"/></td> <td><input type="text" value="3.0"/></td> </tr> <tr> <td>Point 3:</td> <td><input type="text" value="#"/></td> <td><input type="text"/></td> <td><input type="text" value="3.00"/></td> <td><input type="text" value="-2.0"/></td> <td><input type="text" value="3.0"/></td> </tr> <tr> <td>Point 4:</td> <td><input type="text" value="#"/></td> <td><input type="text"/></td> <td><input type="text" value="6.00"/></td> <td><input type="text" value="-2.0"/></td> <td><input type="text" value="3.0"/></td> </tr> <tr> <td>Point 5:</td> <td><input type="text" value="#"/></td> <td><input type="text"/></td> <td><input type="text" value="12.00"/></td> <td><input type="text" value="-2.0"/></td> <td><input type="text" value="3.0"/></td> </tr> <tr> <td>Point 6:</td> <td><input type="text" value="#"/></td> <td><input type="text"/></td> <td><input type="text" value="20.00"/></td> <td><input type="text" value="-2.0"/></td> <td><input type="text" value="3.0"/></td> </tr> <tr> <td>Point 7:</td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>Point 8:</td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>Point 9:</td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>Point 10:</td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td colspan="2">&lt;Point Cal. For Master Curve&gt;</td> <td colspan="2">No. of Correction Points <input type="text"/></td> <td colspan="2">Use Master Curve <input type="text"/></td> </tr> <tr> <td colspan="2"></td> <td colspan="2"></td> <td colspan="2" style="text-align: center;">OD Range</td> </tr> <tr> <td></td> <td>Calibrator</td> <td>OD</td> <td>Conc</td> <td>Low</td> <td>High</td> </tr> <tr> <td>Point-1</td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td>Point-2</td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> <td><input type="text"/></td> </tr> <tr> <td colspan="2">MB Type Factor: <input type="text"/></td> <td colspan="2">1-Point Calibration Point</td> <td colspan="2"><input type="text" value="None"/></td> </tr> </table>						Calibration Type: <input type="text" value="6AB"/>		Formula: <input type="text" value="EIA Type 1"/>		Counts: <input type="text" value="2"/>		<Calibrator Parameters>							Calibrator	OD	Conc	Range						Low	High	Point 1:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="0.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Point 2:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="1.50"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Point 3:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="3.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Point 4:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="6.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Point 5:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="12.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Point 6:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="20.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Point 7:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	Point 8:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	Point 9:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	Point 10:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<Point Cal. For Master Curve>		No. of Correction Points <input type="text"/>		Use Master Curve <input type="text"/>						OD Range			Calibrator	OD	Conc	Low	High	Point-1	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	Point-2	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	MB Type Factor: <input type="text"/>		1-Point Calibration Point		<input type="text" value="None"/>	
Calibration Type: <input type="text" value="6AB"/>		Formula: <input type="text" value="EIA Type 1"/>		Counts: <input type="text" value="2"/>																																																																																																																									
<Calibrator Parameters>																																																																																																																													
	Calibrator	OD	Conc	Range																																																																																																																									
				Low	High																																																																																																																								
Point 1:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="0.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>																																																																																																																								
Point 2:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="1.50"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>																																																																																																																								
Point 3:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="3.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>																																																																																																																								
Point 4:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="6.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>																																																																																																																								
Point 5:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="12.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>																																																																																																																								
Point 6:	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="20.00"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>																																																																																																																								
Point 7:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>																																																																																																																								
Point 8:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>																																																																																																																								
Point 9:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>																																																																																																																								
Point 10:	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>																																																																																																																								
<Point Cal. For Master Curve>		No. of Correction Points <input type="text"/>		Use Master Curve <input type="text"/>																																																																																																																									
				OD Range																																																																																																																									
	Calibrator	OD	Conc	Low	High																																																																																																																								
Point-1	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>																																																																																																																								
Point-2	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>	<input type="text"/>																																																																																																																								
MB Type Factor: <input type="text"/>		1-Point Calibration Point		<input type="text" value="None"/>																																																																																																																									
<table style="width:100%;"> <tr> <td>Slope Check</td> <td><input type="text" value="-"/></td> </tr> <tr> <td>Allowance Range Check</td> <td></td> </tr> <tr> <td><input type="checkbox"/> Reagent Blank</td> <td><input type="text"/></td> </tr> <tr> <td><input type="checkbox"/> Calibration</td> <td><input type="text"/></td> </tr> <tr> <td>Advanced Calibration Operation</td> <td><input type="text" value="No"/></td> </tr> <tr> <td>Interval (RB/ACAL)</td> <td><input type="text"/></td> </tr> <tr> <td><input type="checkbox"/> Lot Calibration</td> <td></td> </tr> </table>						Slope Check	<input type="text" value="-"/>	Allowance Range Check		<input type="checkbox"/> Reagent Blank	<input type="text"/>	<input type="checkbox"/> Calibration	<input type="text"/>	Advanced Calibration Operation	<input type="text" value="No"/>	Interval (RB/ACAL)	<input type="text"/>	<input type="checkbox"/> Lot Calibration																																																																																																											
Slope Check	<input type="text" value="-"/>																																																																																																																												
Allowance Range Check																																																																																																																													
<input type="checkbox"/> Reagent Blank	<input type="text"/>																																																																																																																												
<input type="checkbox"/> Calibration	<input type="text"/>																																																																																																																												
Advanced Calibration Operation	<input type="text" value="No"/>																																																																																																																												
Interval (RB/ACAL)	<input type="text"/>																																																																																																																												
<input type="checkbox"/> Lot Calibration																																																																																																																													
<table style="width:100%;"> <tr> <td>Stability</td> <td>Reagent Blank</td> <td><input type="text" value="#"/></td> <td>Day</td> <td><input type="text" value="#"/></td> <td>Hour</td> </tr> <tr> <td></td> <td>Calibration</td> <td><input type="text" value="#"/></td> <td>Day</td> <td><input type="text" value="#"/></td> <td>Hour</td> </tr> </table>						Stability	Reagent Blank	<input type="text" value="#"/>	Day	<input type="text" value="#"/>	Hour		Calibration	<input type="text" value="#"/>	Day	<input type="text" value="#"/>	Hour																																																																																																												
Stability	Reagent Blank	<input type="text" value="#"/>	Day	<input type="text" value="#"/>	Hour																																																																																																																								
	Calibration	<input type="text" value="#"/>	Day	<input type="text" value="#"/>	Hour																																																																																																																								
<input type="checkbox"/> with Conc-0																																																																																																																													

# User defined

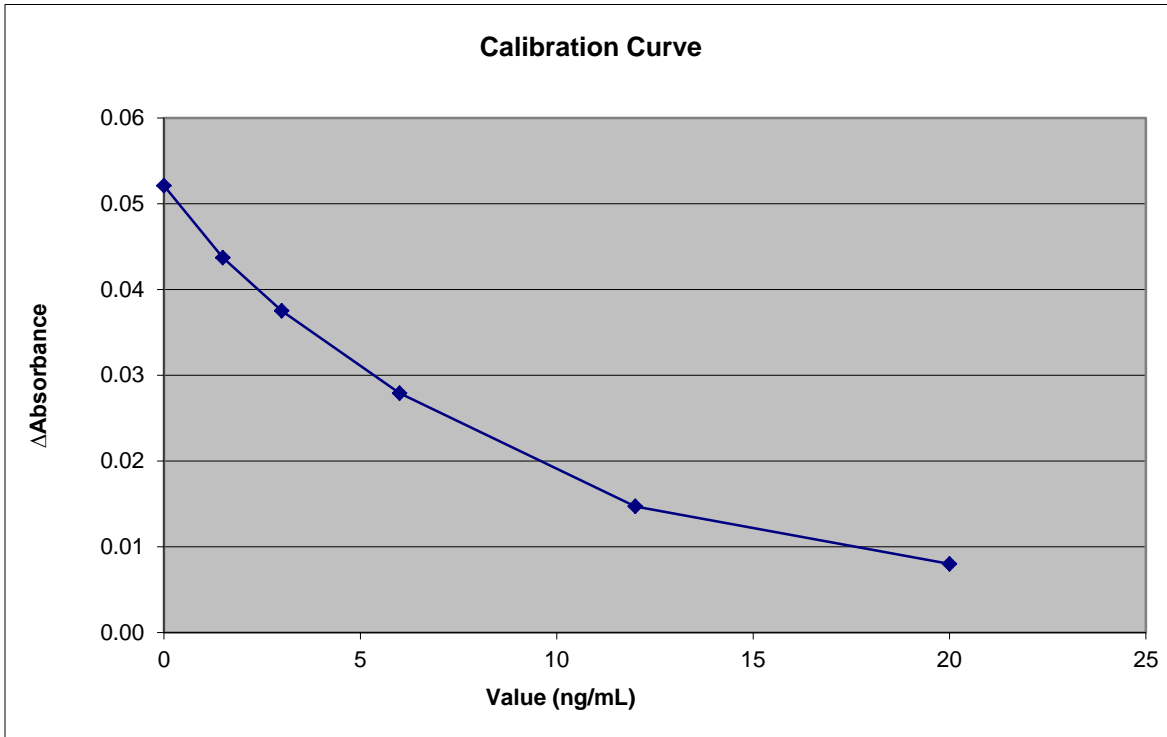


## Results and Data Interpretation

**Performance Data**

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

**Example Calibration Curve, Everolimus (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:

<b>Controls</b>	<b>Control 1</b>	<b>Control 2</b>	<b>Control 3</b>
<b>AU480</b>			
Mean (ng/mL)	3.9	7.5	14.1
Within-Run SD (ng/mL)	0.28	0.46	0.81
Within-Run CV (%)	7.1	6.1	5.7
Total SD (ng/mL)	0.41	0.52	1.02
Total CV (%)	10.5	6.9	7.2
<b>AU680</b>			
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
<b>AU5800</b>			
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

*Continued on next page*

**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 1.5 to 20.0 ng/mL. Error flags will appear for samples recovering above or below the assay range.

The Everolimus assay recovered between 93 – 103% of expected values on the AU480.

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.

**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 AU480 was 0.3 ng/mL.

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 twenty-one blood samples were assayed with the QMS Everolimus Assay on the Beckman Coulter AU480 and tested with reference method Hitachi 917.

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 AU480 =  $0.95 \times (\text{Hitachi 917}) - 0.10$  with a correlation coefficient of 0.992

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.

---

© 2016 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.



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

---

*End*