

**DRI<sup>®</sup> BARBITURATE SERUM TOX APPLICATION**  
**Beckman Coulter AU400<sup>®</sup>, AU480<sup>®</sup>, AU640<sup>®</sup>, AU680<sup>®</sup>,**  
**AU2700<sup>®</sup>, AU5400<sup>®</sup>, AU5800<sup>®</sup>**



Catalog No. 0911

The DRI Barbiturate Serum Tox Assay is intended for the qualitative and semiquantitative determination of barbiturates in human serum or plasma.

For In Vitro Diagnostic Use Only

**Intended Use** 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** Materials available from Microgenics, a part of Thermo Fisher Scientific:

Item	Size	Catalog Number
DRI Barbiturate Serum Tox Kit	R1 25 mL R2 8 mL	0911
DRI Serum Tox Negative Calibrator	1 x 10 mL	0962
DRI Serum Tox Calibrator 1	1 x 5 mL	0963
DRI Serum Tox Calibrator 2	1 x 5 mL	0965
DRI Serum Tox Calibrator 3	1 x 5 mL	0967
DRI Serum Tox Calibrator 4	1 x 5 mL	0976
MAS TOX Control 1 – 3	6 x 5 mL	10011608

To place an order or for technical service, contact:

In USA	In Europe
Tel: (800) 232-3342 Fax: (510) 979-5420	Tel: +49 (0)851-88 6890 Fax: +49 (0)851-88 68910

**Microgenics Corporation, part of Thermo Fisher Scientific**  
**46500 Kato Road, Fremont, CA 94538 USA**  
**U.S. Toll free: (800) 232-3342 / Tel: (510) 979-5000**  
**U.S. Toll free fax: (888) 527-8001 / Fax: (510) 979-5420**

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

*Continued on next page*

---

**Reagent  
Storage**

Refer to the package insert for information on reagent storage.

---

**Analyzer  
Procedure**

Refer to the operator's manuals for information on analyzer operation.

---

**Results and  
Data  
Interpretation**

---

Results for samples will be printed in ng/mL.

# DRI Barbiturate Serum Tox Qualitative Application

## Beckman Coulter System Parameters, AU400/AU640/AU2700/AU5400

Specific Test Parameters	
General	LIH ISE Range
Test Name:	SBARBQ < > Type: Serum Operation: Yes
Sample:	Volume: 4.0 $\mu$ L Dilution: 0 $\mu$ L Pre-Dilution Rate: 1
Reagents:	R1 Volume: 180 $\mu$ L Dilution: 0 $\mu$ L Min OD Max OD
	R2 Volume: 60 $\mu$ L Dilution: 0 $\mu$ L L: -2.00 H: 2.50
Wavelength:	Pri: 340 Sec: 520 Reagent OD limit: First L: -2.00 First H: 2.50
Method:	FIXED1* Last L: -2.00 Last H: 2.50
Reaction slope:	+
Measuring Point 1:	First: 14 Last: 20 Dynamic Range: L: # H: #
Measuring Point 2:	First: Last: Correlation Factor: A: 1 B: 0
Linearity:	%
No-Lag-Time:	No On-board stability period: #

Specific Test Parameters	
General	LIH ISE Range
Test Name:	SBARBQ < > Type: Serum
Value/Flag:	# Level L: # Level H: †
Normal Ranges:	Age L Age H
	Sex Year Month Year Month L H
<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.	None Selected # #
<input type="checkbox"/> 8.	Out of Range L H # #
Panic Value:	# # Unit: % Decimal places: #

Calibration Specific	
General	ISE
Test Name:	SBARBQ < > Type: Serum
Calibration Type:	AA Formula: Y=AX+B Counts: 2 Process: CONC
Point 1:	Cal. No.: # OD: CONC: 0 Factor/OD-L: -99999 Factor/OD-H: 99999
Point 2:	Cal. No.: # OD: CONC: 100 Factor/OD-L: -99999 Factor/OD-H: 99999
Point 3:	Cal. No.: OD: CONC: Factor/OD-L: Factor/OD-H:
Point 4:	Cal. No.: OD: CONC: Factor/OD-L: Factor/OD-H:
Point 5:	Cal. No.: OD: CONC: Factor/OD-L: Factor/OD-H:
Point 6:	Cal. No.: OD: CONC: Factor/OD-L: Factor/OD-H:
Point 7:	Cal. No.: OD: CONC: Factor/OD-L: Factor/OD-H:
1-Point Cal. Point:	<input type="checkbox"/> With CONC-0 Slope Check + Advanced Calibration: #
MB Type Factor:	Calibration Stability Period: #

# User defined  
 \* Can also be run as RATE1

# DRI Barbiturate Serum Tox Semiquantitative Application

## Beckman Coulter System Parameters, AU400/AU640/AU2700/AU5400

Specific Test Parameters											
General		LIH	ISE	Range							
Test Name:		SBARSQ		<	>	Type:	Serum		Operation:	Yes	
Sample:	Volume	4.0	μL	Dilution	0	μL	Pre-Dilution Rate:	1			
Reagents:	R1 Volume	180	μL	Dilution	0	μL	Min OD	Max OD			
	R2 Volume	60	μL	Dilution	0	μL	L	-2.00	H	2.50	
Wavelength:	Pri.	340	▽	Sec.	520	▽	Reagent OD limit:	First L	-2.00	First H	2.50
Method:	FIXED1*					▽	Last L	-2.00	Last H	2.50	
Reaction slope:	+					▽	Dynamic Range:	L	0	H	6000
Measuring Point 1:	First	14	Last	20							
Measuring Point 2:	First		Last								
Linearity:						%	A	1.0000	B	0.0000	
No Lag Time:	No					▽	On-board stability period:	#			

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:		SBARSQ		<	>	Type:	Serum			
Value/Flag:	#		▽	Level L:	#		Level H:	#		
Normal Ranges:		Age L			Age H			L		H
	Sex	Year	Month	Year	Month					
<input type="radio"/>	1.	#	▽	#	#	#	#	#	#	
<input type="radio"/>	2.	#	▽	#	#	#	#	#	#	
<input type="radio"/>	3.	#	▽	#	#	#	#	#	#	
<input type="radio"/>	4.	#	▽	#	#	#	#	#	#	
<input type="radio"/>	5.	#	▽	#	#	#	#	#	#	
<input type="radio"/>	6.	#	▽	#	#	#	#	#	#	
<input type="radio"/>	7.	None Selected					#	#	#	#
<input type="radio"/>	8.	Out of Range		L	H	#	#	#	#	
Panic Value:	#		L	#	H	#	Unit:	ng/mL	Decimal places:	#

Calibration Specific										
General		ISE								
Test Name:		SBARSQ		<	>	Type:	Serum			
Calibration Type:		5AB		▽	Formula:	EIA TYPE 1		▽	Counts:	2
							Process:	CONC		▽
Point 1:	Cal. No.	#	OD		CONC	0	Factor/OD-L	-2.0000	Factor/OD-H	2.5000
Point 2:	Cal. No.	#	OD		CONC	500	Factor/OD-L	-2.0000	Factor/OD-H	2.5000
Point 3:	Cal. No.	#	OD		CONC	1000	Factor/OD-L	-2.0000	Factor/OD-H	2.5000
Point 4:	Cal. No.	#	OD		CONC	3000	Factor/OD-L	-2.0000	Factor/OD-H	2.5000
Point 5:	Cal. No.	#	OD		CONC	6000	Factor/OD-L	-2.0000	Factor/OD-H	2.5000
Point 6:	Cal. No.		OD		CONC		Factor/OD-L		Factor/OD-H	
Point 7:	Cal. No.		OD		CONC		Factor/OD-L		Factor/OD-H	
1-Point Cal. Point:		<input type="checkbox"/>	With CONC-0	Slope Check	+	▽	Advanced Calibration:			▽
MB Type Factor:							Calibration Stability Period:	#		

# User defined  
 \* Can also be run as RATE1

# DRI Barbiturate Serum Tox Qualitative Application

## Beckman Coulter System Parameters, AU480/AU680

Specific Test Parameters

General LIH ISE Range

Test Name: SBARBQ < > Type: Serum Operation: Yes

Sample Volume: 3.6 μL Dilution: 0 μL OD Limit: Min. OD: -2.00 Max. OD: 3.00

Pre-Dilution Rate: 1

Reagents Volume: R1(R1-1) 162 μL Dilution: 0 μL Reagent OD limit: First Low: -2.00 High: 3.00 Last Low: -2.00 High: 3.00

R2 (R2-1) 54 μL Dilution: 0 μL Dynamic Range Low: # High: #

Correlation Factor A: 1 B: 0

Factor for Maker A: 1 B: 0

Wavelength: Pri. 340 nm Sec. 520 nm

Method: FIXED1\*

Reaction slope: + Onboard Stability: 32 Days 0 Hour

Measuring Point 1: First 14 Last 20 LIH Influence Check: #

Measuring Point 2: First Last

Linearity: % Lipemia: Icterus: Hemolysis:

No Lag Time: No

Specific Test Parameters

General ISE Range

Test Name: SBARBQ < > Type: Serum

Value/Flag: # Level L: # Level H: †

Specific Ranges:

	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: % Decimal Places: #

Panic Value: Low # High #

Calibration Specific

General ISE

Test Name: SBARBQ < > Type: Urine  Use Serum Cal.

Calibration Type: AA Formula: Y=AX+B Counts: 2

<Calibrator Parameters>

Calibrator †	OD	Conc	Factor Range	
			Low	High
Point 1: #		0	-99999	999999
Point 2: #		100	-99999	999999
Point 3:				
Point 4:				
Point 5:				
Point 6:				
Point 7:				
Point 8:				
Point 9:				
Point10:				

Slope Check: +

Allowable Range Check:  Reagent Blank  Calibration

Advanced Calibration: Operation Interval (RB/ACAL):

<Point Cal. For Master Curve> No. of Correction Points: Use Master Curve:  Lot Calibration

Calibrator	OD	Conc	OD Range		Stability
			Low	High	Reagent Blanks Calibration
Point 1:					# Day # Hour
Point 2:					18 Day 0 Hour

MB Type Factor: 1-Point Calibration Point:  With CONC-0

# User defined.  
\* Can also be run as RATE1

# DRI Barbiturate Serum Tox Semiquantitative Application

## Beckman Coulter System Parameters, AU480/AU680

Specific Test Parameters											
General		LIH	ISE	Range							
Test Name:	SBARSQ ▾			<	>	Type:	Serum ▾		Operation:	Yes ▾	
Sample Volume	3.6 μL		Dilution	0 μL		OD Limit					
Pre-Dilution Rate	1		Min. OD	-2.000		Max. OD	3.00				
Reagents Volume:	R1(R1-1) 162 μL		Dilution	0 μL		Reagent OD limit:					
			First Low	-2.000		High	3.000				
			Last Low	-2.000		High	3.000				
	R2 (R2-1) 54 μL		Dilution	0 μL		Dynamic Range Low	0		High	6000	
			Correlation Factor A	1		B	0				
Wavelength:	Pri.	340 nm		Sec.	520 nm		Factor for Maker A	1		B	0
Method:	FIXED1* ▾										
Reaction slope:	+ ▾		Onboard Stability	32 Days		0 Hour					
Measuring Point 1:	First	14		Last	20		LIH Influence Check	▾			
Measuring Point 2:	First			Last			Lipemia	▾			
Linearity:	▾ %										
No Lag Time:	No ▾										
			Icterus	▾				▾			
			Hemolysis	▾				▾			

Specific Test Parameters										
General		ISE	Range							
Test Name:	SBARSQ ▾			<	>	Type:	Serum ▾			
Value/Flag:	# ▾		Level L:	# ▾		Level H:	# ▾			
Specific Ranges:										
	From				To			Panic Value		
	Sex	Year	Month	Year	Month	Low	High		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	# ▾						

Calibration Specific										
General		ISE								
Test Name:	SBARSQ ▾			<	>	Type:	Serum ▾		<input type="checkbox"/> Use Serum Cal.	
Calibration Type:	5AB ▾		Formula:	POLYGONAL ▾		Counts:	2			
<Calibrator Parameters>										
Calibrator ↑	OD	Conc	Factor Range				Slope Check			
			Low	High			+ ▾			
Point 1:	# ▾	0	-2.0000	3.0000			Allowable Range Check			
Point 2:	# ▾	500	-2.0000	3.0000			<input type="checkbox"/> Reagent Blank			
Point 3:	# ▾	1000	-2.0000	3.0000			<input type="checkbox"/> Calibration			
Point 4:	# ▾	3000	-2.0000	3.0000			Advanced Calibration			
Point 5:	# ▾	6000	-2.0000	3.0000			Operation			
Point 6:	▾						Interval (RB/ACAL)			
Point 7:	▾						▾			
Point 8:	▾									
Point 9:	▾									
Point10:	▾									
<Point Cal. For Master Curve>										
Calibrator	OD	Conc	No. of Correction Points		Use Master Curve		<input type="checkbox"/> Lot Calibration			
			▾		▾					
			OD Range				Stability			
			Low	High			# Day # Hour			
Point 1:							Reagent Blanks			
Point 2:							Calibration			
							18 Day 0 Hour			
MB Type Factor:			1-Point Calibration Point		▾		<input type="checkbox"/> With CONC-0			

# User Defined  
 \* Can also be run as Rate1

# DRI Barbiturate Serum Tox Qualitative Application

## Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: <input type="text" value="SBARBQ"/> < > Type: <input type="text" value="Serum"/> Operation <input type="text" value="Yes"/>					
Sample Volume	<input type="text" value="3.6"/> $\mu\text{L}$	Dilution	<input type="text" value="0"/> $\mu\text{L}$	OD Limit	
Pre-Dilution Rate	<input type="text" value="1"/>	Diluent Bottle	<input type="text" value=""/>	Min.OD	<input type="text" value="-2.00"/> Max.OD <input type="text" value="3.00"/>
Rgt. Volume	R1(R1-1) <input type="text" value="162"/> $\mu\text{L}$	Dilution	<input type="text" value="0"/> $\mu\text{L}$	Reagent OD Limit	
	R1-2 <input type="text" value=""/>	Dilution	<input type="text" value=""/>	1 <sup>st</sup> . Low	<input type="text" value="-2.00"/> High <input type="text" value="3.00"/>
	R2(R2-1) <input type="text" value="54"/> $\mu\text{L}$	Dilution	<input type="text" value="0"/> $\mu\text{L}$	Last Low	<input type="text" value="-2.00"/> High <input type="text" value="3.00"/>
Common Rgt. Type	<input type="text" value="None"/>	Name	<input type="text" value=""/>	Dynamic Range Low	<input type="text" value="#"/> High <input type="text" value="#"/>
Wavelength	Pri <input type="text" value="340"/> $\text{nm}$	Sec.	<input type="text" value="520"/> $\text{nm}$	Correlation Factor A	<input type="text" value="1"/>
Method	<input type="text" value="FIXED1*"/>			Factor for Maker A	<input type="text" value="1"/>
Reaction Slope	<input type="text" value="+"/>			Onboard Stability Period	<input type="text" value="32"/> Day <input type="text" value="0"/> Hour
Measuring Point1 1 <sup>st</sup>	<input type="text" value="14"/>	Last	<input type="text" value="20"/>	LIH Influence Check	<input type="text" value=""/>
Measuring Point2 1 <sup>st</sup>	<input type="text" value=""/>	Last	<input type="text" value=""/>	Lipemia	<input type="text" value=""/>
Linearity Limit	<input type="text" value=""/>			Icterus	<input type="text" value=""/>
Lag Time Check	<input type="text" value="No"/>			Hemolysis	<input type="text" value=""/>

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: <input type="text" value="SBARBQ"/> < > Type: <input type="text" value="Serum"/>					
Value/Flag:	<input type="text" value="#"/>	Level	Low <input type="text" value="#"/>	High	<input type="text" value="†"/>
Specific Ranges:	From	To	Low	High	
	Sex	Year	Month	Year	Month
o 1.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
o 2.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
o 3.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
o 4.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
o 5.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
o 6.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
7.	Standard demographics				<input type="text" value="#"/>
8.	Not within expected values				<input type="text" value="#"/>
Panic Value	Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>	Unit <input type="text" value="%"/> Decimal Places <input type="text" value="#"/>

Parameters		Calibration Parameters			
Calibrators	Calibration Specific				
General	ISE				
Test Name: <input type="text" value="SBARBQ"/> < > Type <input type="text" value="Serum"/> Cuvette . <input type="text" value=""/>					
<input type="checkbox"/> Use Serum Cal.					
Calibration Type:	<input type="text" value="AA"/>	Formula:	<input type="text" value="Y=AX+B"/>		Counts: <input type="text" value="2"/>
<Calibrator Parameters>					
	Calibrator	OD	Conc	Range	Slope Check <input type="text" value="+"/>
Point 1:	<input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="0"/>	Low <input type="text" value="-99999"/> High <input type="text" value="999999"/>	Allowance Range Check
Point 2:	<input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="100"/>	Low <input type="text" value="-99999"/> High <input type="text" value="999999"/>	<input type="checkbox"/> Reagent Blank <input type="text" value=""/>
Point 3:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		<input type="checkbox"/> Calibration <input type="text" value=""/>
Point 4:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		Advanced Calibration
Point 5:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		Operation <input type="text" value="No"/>
Point 6:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		Interval (RB/ACAL) <input type="text" value=""/>
Point 7:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		
Point 8:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		
Point 9:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		
Point 10:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>		
<Point Cal. For	No. of Correction Points	<input type="text" value=""/>	Use Master Curve	<input type="text" value=""/>	<input type="checkbox"/> Lot Calibration
Master Curve>			OD Range		
	Calibrator	OD	Conc	Low	High
Point-1	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>
Point-2	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>
MB Type Factor:	<input type="text" value=""/>	1-Point Calibration Point	<input type="text" value="None"/>	<input type="checkbox"/> with Conc-0	
				Stability	Reagent Blank <input type="text" value="#"/> Day <input type="text" value="#"/> Hour
					Calibration <input type="text" value="18"/> Day <input type="text" value="0"/> Hour

# User defined.  
\* Can also be run as RATE1

# DRI Barbiturate Serum Tox Semiquantitative Application

## Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: SBARSQ < > Type: Serum Operation Yes					
Sample Volume	3.6	μL	Dilution	0	μL
Pre-Dilution Rate	1	∇	Diluent Bottle	#	∇
Rgt. Volume	R1(R1-1)	162	μL	Dilution	0
	R1-2		μL	Dilution	
	R2(R2-1)	54	μL	Dilution	0
Common Rgt. Type	None	Name	None	Correlation Factor A	1
Wavelength	Pri 340	∇nm	Sec. 520	∇nm	Factor for Maker A 1
Method	FIXED1*	∇			
Reaction Slope	+	∇			
Measuring Point1	1 <sup>st</sup>	14			
Measuring Point2	1 <sup>st</sup>				
Linearity Limit		%			
Lag Time Check	No	∇			
				Dynamic Range Low	0
				High	1000
				Onboard Stability Period	32 Day 0 Hour
				LIH Influence Check	# ∇
				Lipemia	∇
				Icterus	∇
				Hemolysis	∇

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: SBARSQ < > Type: Serum					
Value/Flag:	#	∇			
Specific Ranges:	From	Level	To	Low	High
	Sex	Year	Month	Year	Month
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 #

Parameters		Calibration Parameters			
Calibrators	Calibration Specific				
General	ISE				
Test Name: SBARSQ < > Type: Serum Cuvette .					
<input type="checkbox"/> Use Serum Cal.					
Calibration Type:	5AB	∇	Formula:	POLYGONAL	∇
			Counts:	2	∇
<Calibrator Parameters>					
Calibrator	OD	Conc	Low	High	Slope Check
Point 1:	#	0	-2.00	3.00	+
Point 2:	#	500	-2.00	3.00	
Point 3:	#	1000	-2.00	3.00	
Point 4:	#	3000	-2.00	3.00	
Point 5:	#	6000	-2.00	3.00	
Point 6:					
Point 7:					
Point 8:					
Point 9:					
Point 10:					
Allowance Range Check					
<input type="checkbox"/> Reagent Blank					
<input type="checkbox"/> Calibration					
Advanced Calibration Operation					
Interval (RB/ACAL)					
<Point Cal. For Master Curve>					
No. of Correction Points			Use Master Curve		<input type="checkbox"/> Lot Calibration
Calibrator	OD	Conc	Low	High	Stability
Point-1					Reagent Blank # Day # Hour
Point-2					Calibration 18 Day 0 Hour
MB Type Factor:	1-Point Calibration Point None ∇ <input type="checkbox"/> with Conc-0				

# User defined.  
\* Can also be run as RATE1



**Precision**

Tests for within-run and total precision, evaluated with packaged reagents, controls and calibrators, yielded the following results (N=80/level):

Controls	Control 1	Control 2	Control 3	Cutoff
<b>Qualitative AU480</b>				
Mean (mA/min)	565	639	702	622
Within-Run SD	1.5	2.1	1.7	1.5
Within-Run CV	0.3%	0.3%	0.2%	0.2%
Total SD (A/min)	2.4	2.8	3.1	2.3
Total CV (%)	0.4%	0.4%	0.4%	0.4%
<b>Qualitative AU640*</b>				
Mean (mA/min)	614	N/A	743	656
Within-Run	3.6	N/A	4.8	4.2
Within-Run CV	0.6%	N/A	0.7	0.6%
Total SD (mA/min)	4.3	N/A	4.8	4.2
Total CV (%)	0.7%	N/A	0.7	0.6%
<b>Qualitative AU680</b>				
Mean (mA/min)	1026	1158	1270	1130
Within-Run SD	5.1	6.7	8.4	5.0
Within-Run CV	0.5%	0.6%	0.7%	0.4%
Total SD (mA/min)	7.3	8.4	9.1	7.5
Total CV (%)	0.7%	0.7%	0.7%	0.7%
<b>Qualitative AU5800</b>				
Mean (mA/min)	553	626	690	612
Within-Run SD	8.3	7.4	7.9	8.1
Within-Run CV	1.5%	1.2%	1.1%	1.3%
Total SD (mA/min)	8.8	8.3	9.3	8.9
Total CV (%)	1.6%	1.3%	1.3%	1.5%

Controls	Control 1	Control 2	Control 3	Cutoff
<b>Semiquantitative AU480</b>				
Mean (ng/mL)	362	1399	3340	962
Within-Run SD	8.9	58.8	102.1	22.7
Within-Run CV	2.4%	4.2%	3.1%	2.4%
Total SD (ng/mL)	16.5	76.6	182.1	33.4
Total CV (%)	4.6%	5.5%	5.5%	3.5%
<b>Semiquantitative AU640*</b>				
Mean (ng/mL)	579	N/A	2961	1032
Within-Run SD	35.4	N/A	135.3	71.5
Within-Run CV	6.1%	N/A	4.6%	6.9%
Total SD (ng/mL)	42.6	N/A	140.6	72.6
Total CV (%)	7.4%	N/A	4.8%	7.0%
<b>Semiquantitative AU680</b>				
Mean (ng/mL)	348	1424	3256	1026
Within-Run SD	18.0	97.8	192.7	57.9
Within-Run CV	5.2%	6.9%	5.9%	5.6%
Total SD (ng/mL)	26.4	136.2	226.4	94.9
Total CV (%)	7.6%	9.6%	7.0%	9.3%
<b>Semiquantitative AU5800</b>				
Mean (ng/mL)	354	1417	3278	1060
Within-Run SD	24.0	97.5	208.4	74.3
Within-Run CV	6.8%	6.9%	6.4%	7.0%
Total SD (ng/mL)	24.3	113.2	245.5	88.2
Total CV (%)	6.9%	8.0%	7.5%	8.3%

AU640\* N=60/level

**Accuracy and Correlation**

Patient samples were assayed using the DRI Barbiturate Serum Tox Assay on the Beckman Coulter AU480, AU640, AU680, and AU5800; and tested against the reference analyzer, the Hitachi 717.

Qualitative Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	94.8	90.5	93.0
AU640	100.0	98.7	99.0
AU680	100.0	97.9	99.0
AU5800	97.1	90.7	94.7

**DRI Serum Tox Barbiturate Assay Qualitative Method Comparison**

		Hitachi 717				Hitachi 717	
		+	-			+	-
AU480	+	55	4	AU640	+	25	1
	-	3	38		-	0	74
AU680	+	52	1	AU5800	+	68	4
	-	0	47		-	2	39

Semiquantitative Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	100.0	95.5	98.0
AU640	96.3	100.0	99.0
AU680	100.0	95.5	98.0
AU5800	100.0	90.9	96.0

**DRI Serum Tox Barbiturate Assay Semiquantitative Method Comparison**

		Hitachi 717				Hitachi 717	
		+	-			+	-
AU480	+	56	2	AU640	+	26	0
	-	0	42		-	1	73
AU680	+	56	2	AU5800	+	56	4
	-	0	42		-	0	40

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

End