

DRITM OXYCODONE APPLICATION
BECKMAN COULTER AU480[®], AU680[®], AU5800[®]



Catalog No. 100248, 100249, 10015632

The DRI Oxycodone Enzyme Immunoassay is intended for the qualitative and semiquantitative determination of oxycodone in human urine.



For In Vitro Diagnostic Use Only
Rx 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. For package inserts, visit www.thermoscientific.com/diagnostics and enter the assay name in the *Search* field.

Ordering Information

For ordering information, visit www.thermoscientific.com/diagnostics and enter the assay name in the *Search* field. Not all intended uses and applications mentioned here are available in every country. Please contact your local sales representative or distributor for more information.

For Technical Service, contact:

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



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.

Dispense adequate amounts of Antibody/Substrate Reagent A (first reagent) and Enzyme Conjugate Reagent E (second reagent) into appropriate containers. **Make sure that the reagents have equilibrated to the temperature of the analyzer reagent compartment before starting analysis.**

Calibration

Refer to package insert for information on calibration.

**Results and
Data
Interpretation**

Refer to package insert for information on results and data interpretation.

DRI Oxycodone Assay – Qualitative (100 ng/mL cutoff) Beckman Coulter System Parameters, AU480 / AU680

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾	Operation	Yes ▾		
Sample Volume	13.9 μL	Dilution	0 ▾ μL	OD Limit		Min.OD	-2.0000	Max.OD	3.0000
Pre-Dilution Rate	1 ▾								
Reagent Volume	R1(R1-1) 87 μL	Dilution	0 μL	Reagent OD Limit		First	Low -2.0000	High	3.0000
	R2(R2-1) 87 μL	Dilution	0 μL			Last	Low -2.0000	High	3.0000
Common Reagent	Type	None	Name		Dynamic Range	Low	#	High	#
Wavelength	Pri.	340 ▾ nm	Sec.	410 ▾ nm	Correlation Factor	A	1	B	†
Method		FIXED* ▾			Factor for Maker	A	1	B	0
Reaction slope		+ ▾							
Measuring Point-1	First	17	Last	21	Onboard Stability Period		30 Day	# Hour	
Measuring Point-2	First		Last						
Linearity Limit									
Lag Time Check		No ▾							

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾				
Value/Flag	# ▾								
Level		Low	-9999999	High	‡	Panic Value			
						Low	#	High	#
Specific Ranges:		From		To					
	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	#						

User defined.

* Can also be run as RATE

- Option 1: Run a reagent blank (blue rack). Run the cutoff calibrator in a white rack. Compare the sample response to the cutoff calibrator response to determine if the sample is positive or negative. Positive samples will not be flagged.
- Option 2: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to 100.
- Option 3: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to zero.

DRI Oxycodone Assay – Qualitative (300 ng/mL cutoff) Beckman Coulter System Parameters, AU480/ AU680

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾	Operation	Yes ▾		
Sample Volume	10.4 μL	Dilution	0 ▾ μL	OD Limit		Min.OD	-2.0000	Max.OD	3.0000
Pre-Dilution Rate	1 ▾								
Reagent Volume	R1(R1-1) 87 μL	Dilution	0 μL	Reagent OD Limit		First	Low -2.0000	High	3.0000
	R2(R2-1) 87 μL	Dilution	0 μL			Last	Low -2.0000	High	3.0000
Common Reagent	Type	None	Name		Dynamic Range	Low	#	High	#
Wavelength	Pri.	340 ▾ nm	Sec.	410 ▾ nm	Correlation Factor	A	1	B	†
Method		FIXED* ▾			Factor for Maker	A	1	B	0
Reaction slope		+ ▾							
Measuring Point-1	First	17	Last	21	Onboard Stability Period		30 Day	# Hour	
Measuring Point-2	First		Last						
Linearity Limit									
Lag Time Check		No ▾							

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾				
Value/Flag	# ▾								
Level		Low	-9999999	High	‡	Panic Value			
						Low	#	High	#
Specific Ranges:		From		To					
	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	#						

User defined

* Can also be run as RATE

† Option 1: Enter 0.0 Option 2: Enter 0.0 Option 3: Enter -100

‡ Option 1: Enter 9999999 Option 2: Enter 100 Option 3: Enter 0.0

- Option 1: Run a reagent blank (blue rack). Run the cutoff calibrator in a white rack. Compare the sample response to the cutoff calibrator response to determine if the sample is positive or negative. Positive samples will not be flagged.
- Option 2: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to 100.
- Option 3: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to zero.

DRI Oxycodone Assay – Qualitative (100 or 300 ng/mL cutoff) Beckman Coulter System Parameters, AU480 / AU680, *continued*

(Option 1)

Calibration Specific									
General		ISE							
Test Name:		#	<	>	Type	Urine	<input type="checkbox"/> Use Serum Cal.		
Calibration Type:		MB	Formula:		Y=AX+B	Counts:	2		
<Calibrator Parameters>									
	Calibrator †	OD	Conc	Factor Range		Slope Check			
				Low	High				
Point 1:	▽					Allowable Range Check			
Point 2:	▽					<input type="checkbox"/> Reagent Blank			
Point 3:	▽					<input type="checkbox"/> Calibration			
Point 4:	▽					Advanced Calibration			
Point 5:	▽					Operation		No	
Point 6:	▽					Interval (RB/ACAL)			
Point 7:	▽								
Point 8:	▽								
Point 9:	▽								
Point10:	▽								
<Point Cal. For Master Curve>									
	Calibrator	OD	Conc	OD Range		Stability			
				Low	High	Reagent Blanks Calibration			
Point 1:						#		Day	
Point 2:						#		Hour	
MB Type Factor:		1000		1-Point Calibration Point		<input type="checkbox"/> With CONC-0			

(Option 2 or 3)

Calibration Specific									
General		ISE							
Test Name:		#	<	>	Type	Urine	<input type="checkbox"/> Use Serum Cal.		
Calibration Type:		AB	Formula:		Y=AX+B	Counts:	2		
<Calibrator Parameters>									
	Calibrator †	OD	Conc	Factor Range		Slope Check			
				Low	High				
Point 1:	▽		100	--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		No	
Point 6:	▽					Interval (RB/ACAL)			
Point 7:	▽								
Point 8:	▽								
Point 9:	▽								
Point10:	▽								
<Point Cal. For Master Curve>									
	Calibrator	OD	Conc	OD Range		Stability			
				Low	High	Reagent Blanks Calibration			
Point 1:						#		Day	
Point 2:						#		Hour	
MB Type Factor:				1-Point Calibration Point		<input type="checkbox"/> With CONC-0			

User defined.

DRI Oxycodone Assay – Semiquantitative (100 and 300 ng/mL cutoff) Beckman Coulter System Parameters, AU480 /AU680

Specific Test Parameters												
General		LIH	ISE	Range								
Test Name:		#	<	>	Type:	Urine	Operation:		Yes			
Sample Volume	10.4	μL	Dilution	0	μL	OD Limit						
Pre-Dilution Rate	1				Min. OD	-2.00	Max. OD	3.00				
Reagents Volume:	R1(R1-1)	87	μL	Dilution	0	μL	Reagent OD limit:					
							First Low	-2.00	Hig	3.00		
							Last Low	-2.00	Hig	3.00		
	R2 (R2-1)	87	μL	Dilution	0	μL	Dynamic Range Low		#	Hig	#	
Wavelength:	Pri.	340	nm	Sec.	410	nm	Correlation Factor A		1	B	0	
Method:	FIXED1						Factor for Maker A		1	B	0	
Reaction slope:	+						Onboard Stability		<input type="checkbox"/>	Days	<input type="checkbox"/>	Hour
Measuring Point 1:	First	17	Last	21			LIH Influence Check		#			
Measuring Point 2:	First		Last				Lipemia		<input type="checkbox"/>			
Linearity:							Icterus		<input type="checkbox"/>			
No Lag Time:							Hemolysis		<input type="checkbox"/>			

Specific Test Parameters										
General		ISE	Range							
Test Name:		OXY100SQ	<	>	Type:	Urine				
Value/Flag:	VALUE	Level L:	#	Level H:	#					
Specific Ranges:										
		From	To		Low	High				
<input type="checkbox"/>	1.	Sex	Year	Month	Year	Month	Low	High		
<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			
						#	#			

User defined.

**DRI Oxycodone Assay – Semiquantitative (100 and 300 ng/mL cutoff)
Beckman Coulter System Parameters, AU480 /AU680, *continued***

Calibration Specific										
General		ISE								
Test Name:		#	<	>	Type	Urine	<input type="checkbox"/> Use Serum Cal.			
Calibration Type:		5AB	Formula:		Polygonal*	Counts:		2		
<Calibrator Parameters>					Slope Check					+
	Calibrator †	OD	Conc	Factor Range		Allowable Range Check				
				Low	High	<input type="checkbox"/> Reagent Blank				
Point 1:	#		0.0	-2.00	3.00	<input type="checkbox"/> Calibration				
Point 2:	#		100.0	-2.00	3.00	Advanced Calibration				
Point 3:	#		300.0	-2.00	3.00	Operation				
Point 4:	#		500.0	-2.00	3.00	Interval (RB/ACAL)				
Point 5:	#		1000.0	-2.00	3.00	No				
Point 6:										
Point 7:										
Point 8:										
Point 9:										
Point10:										
<Point Cal. For Master Curve>					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		None	<input type="checkbox"/> With CONC-0				

User Defined

* Can also be run with EIA Type1

DRI Oxycodone Assay – Qualitative (100 ng/mL cutoff) Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾	Operation	Yes ▾		
Sample Volume	13.9 μL	Dilution	0 ▾ μL	OD Limit		Min.OD	-2.0000	Max.OD	3.0000
Pre-Dilution Rate	1 ▾								
Reagent Volume	R1(R1-1) 87 μL	Dilution	0 μL	Reagent OD Limit		First	Low -2.0000	High	3.0000
	R2(R2-1) 87 μL	Dilution	0 μL			Last	Low -2.0000	High	3.0000
Common Reagent	Type	None	Name		Dynamic Range	Low	#	High	#
Wavelength	Pri.	340 ▾ nm	Sec.	410 ▾ nm	Correlation Factor	A	1	B	†
Method		FIXED* ▾			Factor for Maker	A	1	B	0
Reaction slope		+ ▾							
Measuring Point-1	First	17	Last	21	Onboard Stability Period		30 Day	#	Hour
Measuring Point-2	First		Last						
Linearity Limit									
Lag Time Check		No ▾							

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾				
Value/Flag	# ▾								
Level		Low	-9999999	High	‡	Panic Value			
						Low	#	High	#
Specific Ranges:		From		To					
	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	#						

User defined

* Can also be run as RATE

† Option 1: Enter 0.0 Option 2: Enter 0.0 Option 3: Enter -100

‡ Option 1: Enter 9999999 Option 2: Enter 100 Option 3: Enter 0.0

- Option 1: Run a reagent blank (blue rack). Run the cutoff calibrator in a white rack. Compare the sample response to the cutoff calibrator response to determine if the sample is positive or negative. Positive samples will not be flagged.
- Option 2: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to 100.
- Option 3: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to zero.

DRI Oxycodone Assay – Qualitative (300 ng/mL cutoff) Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾	Operation	Yes ▾		
Sample Volume	10.4 μL	Dilution	0 ▾ μL	OD Limit		Min.OD	-2.0000	Max.OD	3.0000
Pre-Dilution Rate	1 ▾								
Reagent Volume	R1(R1-1) 87 μL	Dilution	0 μL	Reagent OD Limit		First	Low -2.0000	High	3.0000
	R2(R2-1) 87 μL	Dilution	0 μL			Last	Low -2.0000	High	3.0000
Common Reagent	Type	None	Name		Dynamic Range	Low	#	High	#
Wavelength	Pri.	340 ▾ nm	Sec.	410 ▾ nm	Correlation Factor	A	1	B	†
Method		FIXED* ▾			Factor for Maker	A	1	B	0
Reaction slope		+ ▾							
Measuring Point-1	First	17	Last	21	Onboard Stability Period		30 Day	#	Hour
Measuring Point-2	First		Last						
Linearity Limit									
Lag Time Check		No ▾							

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾				
Value/Flag	# ▾								
Level		Low	-9999999	High	‡	Panic Value			
						Low	#	High	#
Specific Ranges:		From		To					
	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	#						

User defined

* Can also be run as RATE

† Option 1: Enter 0.0 Option 2: Enter 0.0 Option 3: Enter -100

‡ Option 1: Enter 9999999 Option 2: Enter 100 Option 3: Enter 0.0

- Option 1: Run a reagent blank (blue rack). Run the cutoff calibrator in a white rack. Compare the sample response to the cutoff calibrator response to determine if the sample is positive or negative. Positive samples will not be flagged.
- Option 2: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to 100.
- Option 3: Run a reagent blank (blue rack). Calibrate by placing the appropriate cutoff calibrator in the assigned position in the calibration rack (yellow rack). Positive samples will be flagged (P) and will be greater than or equal to zero.

DRI Oxycodone Assay – Qualitative (100 or 300 ng/mL cutoffs) Beckman Coulter System Parameters, AU5800, *continued*

(Option 1)

Parameters	Calibration Parameters		
Calibrators	Calibration Specific	STAT Table Calibration	
Test Name <input type="text" value="#"/> <input type="button" value="v"/> <input type="button" value="<"/> <input type="button" value=">"/>	Type <input type="text" value="Urine"/> <input type="button" value="v"/>	<input type="checkbox"/> Use Serum Cal.	
Calibration Type <input type="text" value="MB"/> <input type="button" value="v"/>	Formula <input type="text" value="Y=AX+B"/> <input type="button" value="v"/>	Counts <input type="text" value="2"/> <input type="button" value="v"/>	
< Calibrator Parameters >		Range	Slope Check <input type="text" value=""/> <input type="button" value="v"/>
	Calibrator	OD	Conc
			Low High
Point-1	<input type="text" value="v"/>		
Point-2	<input type="text" value="v"/>		
Point-3	<input type="text" value="v"/>		
Point-4	<input type="text" value="v"/>		
Point-5	<input type="text" value="v"/>		
Point-6	<input type="text" value="v"/>		
Point-7	<input type="text" value="v"/>		
Point-8	<input type="text" value="v"/>		
Point-9	<input type="text" value="v"/>		
Point-10	<input type="text" value="v"/>		
< Point Cal. For Master Curve >		No. of Correction Points <input type="text" value="v"/>	Use Master Curve <input type="text" value="v"/>
	Calibrator	OD	Conc
			Low High
Point-1	<input type="text" value="v"/>		
Point-2	<input type="text" value="v"/>		
MB Type Factor <input type="text" value="1000"/>	1-Point Calibration Point <input type="text" value="v"/>	<input type="checkbox"/> with Conc-0	
		Stability	Reagent Blank <input type="text" value=""/> Day <input type="text" value=""/> Hour
		Calibration	<input type="text" value=""/> Day <input type="text" value=""/> Hour
		Advanced Calibration	Operation Interval (RB/ACAL) <input type="text" value="No"/> <input type="button" value="v"/>
			<input type="checkbox"/> Lot Calibration
		Allowable Range Check	<input type="text" value=""/>
		<input type="checkbox"/> Reagent Blank	<input type="text" value=""/>
		<input type="checkbox"/> Calibration	<input type="text" value=""/>

(Option 2 or 3)

Parameters	Calibration Parameters		
Calibrators	Calibration Specific	STAT Table Calibration	
Test Name <input type="text" value="#"/> <input type="button" value="v"/> <input type="button" value="<"/> <input type="button" value=">"/>	Type <input type="text" value="Urine"/> <input type="button" value="v"/>	<input type="checkbox"/> Use Serum Cal.	
Calibration Type <input type="text" value="AB"/> <input type="button" value="v"/>	Formula <input type="text" value="Y=AX+B"/> <input type="button" value="v"/>	Counts <input type="text" value="2"/> <input type="button" value="v"/>	
< Calibrator Parameters >		Factor Range	Slope Check <input type="text" value=""/> <input type="button" value="v"/>
	Calibrator	OD	Conc
			Low High
Point-1	<input type="text" value="#"/> <input type="button" value="v"/>		100 -9999999 9999999
Point-2	<input type="text" value="v"/>		
Point-3	<input type="text" value="v"/>		
Point-4	<input type="text" value="v"/>		
Point-5	<input type="text" value="v"/>		
Point-6	<input type="text" value="v"/>		
Point-7	<input type="text" value="v"/>		
Point-8	<input type="text" value="v"/>		
Point-9	<input type="text" value="v"/>		
Point-10	<input type="text" value="v"/>		
< Point Cal. For Master Curve >		No. of Correction Points <input type="text" value="v"/>	Use Master Curve <input type="text" value="v"/>
	Calibrator	OD	Conc
			Low High
Point-1	<input type="text" value="v"/>		
Point-2	<input type="text" value="v"/>		
MB Type Factor <input type="text" value=""/>	1-Point Calibration Point <input type="text" value="v"/>	<input type="checkbox"/> with Conc-0	
		Stability	Reagent Blank <input type="text" value="#"/> Day <input type="text" value="#"/> Hour
		Calibration	<input type="text" value="#"/> Day <input type="text" value="#"/> Hour
		Advanced Calibration	Operation Interval (RB/ACAL) <input type="text" value="No"/> <input type="button" value="v"/>
			<input type="checkbox"/> Lot Calibration
		Allowable Range Check	<input type="text" value=""/>
		<input type="checkbox"/> Reagent Blank	<input type="text" value=""/>
		<input type="checkbox"/> Calibration	<input type="text" value=""/>

User defined

DRI Oxycodone Assay – Semiquantitative (100 and 300 ng/mL cutoffs) Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾	Operation	Yes ▾		
Sample Volume	10.4 μL	Dilution	0 ▾ μL	OD Limit	Min.OD	-2.0000	Max.OD	3.0000	
Pre-Dilution Rate	1 ▾								
Reagent Volume R1(R1-1)	87 μL	Dilution	0 μL	Reagent OD Limit	First	Low	-2.0000	High	3.0000
					Last	Low	-2.0000	High	3.0000
	R2(R2-1)	87 μL	Dilution	0 μL					
Common Reagent Type	None	Name		Dynamic Range	Low	#	High	#	
Wavelength Pri.	340 ▾ nm	Sec.	410 ▾ nm	Correlation Factor	A	1	B	0	
Method	FIXED1* ▾			Factor for Maker	A	1	B	0	
Reaction slope	+ ▾								
Measuring Point-1	First	17	Last	21	On-board Stability Period	30	Day	#	Hour
Measuring Point-2	First		Last						
Linearity Limit									
Lag Time Check	No ▾								

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	# ▾	<	>	Type	Urine ▾				
Value/Flag Level	# ▾			Low	#	High	#		
						Panic Value			
						Low	#	High	#
Specific Ranges:	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		#	#				
Unit	#	Decimal Places	#						

User defined
* Can also be run as RATE1

**DRI Oxycodone Assay – Semiquantitative (100 and 300 ng/mL cutoffs)
Beckman Coulter System Parameters, AU5800, *continued***

Parameters		Calibration Parameters			
Calibrators		Calibration Specific		STAT Table Calibration	
Test Name	# ▾	<	>	Type	Urine ▾ <input type="checkbox"/> Use Serum Cal.
<div style="border: 1px solid black; padding: 5px;"> Calibration Type 5AB ▾ Formula POLYGONAL* ▾ Counts 2 ▾ </div>					
< Calibrator Parameters >			OD Range		Slope Check
	Calibrator	OD	Conc	Low	High
Point-1	# ▾		0	-2.0000	3.0000
Point-2	# ▾		100	-2.0000	3.0000
Point-3	# ▾		300	-2.0000	3.0000
Point-4	# ▾		500	-2.0000	3.0000
Point-5	# ▾		1000	-2.0000	3.0000
Point-6	▾				
Point-7	▾				
Point-8	▾				
Point-9	▾				
Point-10	▾				
			Slope Check		+ ▾
			Allowable Range Check		
			<input type="checkbox"/> Reagent Blank		
			<input type="checkbox"/> Calibration		
			Advanced Calibration		
			Operation		No ▾
			Interval (RB/ACAL)		▾
			<input type="checkbox"/> Lot Calibration		
< Point Cal. For Master Curve > No. of Correction Points ▾ Use Master Curve ▾					
			OD Range		
	Calibrator	OD	Conc	Low	High
Point-1	▾				
Point-2	▾				
Stability			Reagent Blank		 Day Hour
			Calibration		 Day Hour
MB Type Factor 		1-Point Calibration Point None ▾		<input type="checkbox"/> with Conc-0	

User defined

* Can also be run with EIA type 1.

© 2017 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