

DRI[®] ETHYL GLUCURONIDE APPLICATION **CE**
**BECKMAN COULTER AU400[®], AU480[®], AU640[®], AU680[®],
AU2700[®], AU5400[®], AU5800[®]**

Catalog No. 10011297, 10015626

The Thermo Scientific™ DRI Ethyl Glucuronide Enzyme Immunoassay is intended for the qualitative and semiquantitative determination of ethyl glucuronide (EtG) in human urine at cutoffs of 500 ng/mL and 1000 ng/mL.

For In Vitro Diagnostic Use Only
For Export Only, Not Available in the United States

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 No.
DRI Ethyl Glucuronide Assay	68 mL Kit	10011297
	3 x 18 mL Kit	10015626
DRI Ethyl Glucuronide Negative Calibrator	25 mL	10011207
DRI Ethyl Glucuronide 100 ng/mL Calibrator	10 mL	10011208
DRI Ethyl Glucuronide 500 ng/mL Calibrator	10 mL	10011210
DRI Ethyl Glucuronide 1000 ng/mL Calibrator	10 mL	10011212
DRI Ethyl Glucuronide 2000 ng/mL Calibrator	10 mL	10011213
DRI Ethyl Glucuronide 375 ng/mL Control	25 mL	10012135
DRI Ethyl Glucuronide 625 ng/mL Control	25 mL	10012136
DRI Ethyl Glucuronide 750 ng/mL Control	25 mL	10012137
DRI Ethyl Glucuronide 1250 ng/mL Control	25 mL	10012138

To place an order or for technical service contact:



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 Microgenics GmbH, Spitalhofstrasse 94, D-94032 Passau, Germany

Continued on next page

Reagent Storage

Refer to package insert for information on reagent storage.

Procedure for Analyzer

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

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

Note: Under Specific Test Parameters/General Tab, Linearity % should be left blank, as reflected in the following pages. **Do Not Enter 0.**

Results and Data Interpretation

Results for samples will be printed in ng/mL for the semiquantitative assay and in % units for the qualitative assay.

DRI Ethyl Glucuronide – Qualitative

Beckman Coulter System Parameters, AU400/AU640

Specific Test Parameters			
General	LIH	ISE	Range
Test Name:	ETGQ	<	>
Type:	Urine	Operation:	Yes
Sample:	Volume	35.0	μL Dilution
Reagents:	R1 Volume	80	μL Dilution
	R2 Volume	80	μL Dilution
Wavelength:	Pri.	340	Sec. 410
Method:	RATE*		
Reaction slope:	+		
Measuring Point 1:	First	13	Last 17
Measuring Point 2:	First		Last
Linearity			
No-Lag-Time:	No		
Pre-Dilution Rate:	1		
Min OD	Max OD		
	L	-2.00	H 2.50
Reagent OD limit:	First L	-2.00	First H 2.50
	Last L	-2.00	Last H 2.50
Dynamic Range:	L	#	H #
Correlation Factor:	A	1	B 0
On-board stability period:	#		

Specific Test Parameters			
General	LIH	ISE	Range
Test Name:	ETGQ	<	>
Type:	Urine		
Value/Flag:	#	Level L:	#
		Level H:	#
Normal Ranges:	Age L	Age H	
	Sex	Year	Month
	Year	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.	#	#	#
7.	None Selected		
8.	Out of Range		
Panic Value:	L	H	
	#	#	
Unit:	#	Decimal places:	#

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

User defined
 * Can also be run as FIXED

DRI Ethyl Glucuronide – Semiquantitative Beckman Coulter System Parameters, AU400/AU640

Specific Test Parameters	
General	LIH ISE Range
Test Name:	ETGSQ < > Type: Urine < Operation: Yes <
Sample: Volume	35.0 μL Dilution 0 μL Pre-Dilution Rate: 1
Reagents: R1 Volume	80 μL Dilution 0 μL Min OD Max OD
R2 Volume	80 μL Dilution 0 μL L -2.00 H 2.50
Wavelength: Pri.	340 < Sec. 410 < Reagent OD limit: First L -2.00 First H 2.50
Method:	RATE1* < Last L -2.00 Last H 2.50
Reaction slope:	+ < Dynamic Range: L 0 H 2000
Measuring Point 1:	First 13 Last 17 Correlation Factor: A 1 B 0
Measuring Point 2:	First Last On-board stability period: #
Linearity:	%
No Lag Time:	No <

Specific Test Parameters	
General	LIH ISE Range
Test Name:	ETGSQ < > Type: Urine <
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.	# < # # # # # # #
7.	None Selected
8.	Out of Range L H
Panic Value:	# # Unit: ng/mL Decimal places: #

Calibration Specific	
General	ISE
Test Name:	ETGSQ < > Type: Urine <
Calibration Type:	5AB < Formula: POLYGONAL < Counts: 2 Process: CONC <
Point 1:	Cal. No. 1 OD CONC 0.0 Factor/OD-L -2.0 Factor/OD-H 2.5
Point 2:	2 OD CONC 100.0 Factor/OD-L -2.0 Factor/OD-H 2.5
Point 3:	3 OD CONC 500.0 Factor/OD-L -2.0 Factor/OD-H 2.5
Point 4:	4 OD CONC 1000.0 Factor/OD-L -2.0 Factor/OD-H 2.5
Point 5:	5 OD CONC 2000.0 Factor/OD-L -2.0 Factor/OD-H 2.5
Point 6:	OD CONC Factor/OD-L Factor/OD-H
Point 7:	OD CONC Factor/OD-L Factor/OD-H
1-Point Cal. Point:	With CONC-0
MB Type Factor:	Calibration Stability Period: #

User defined
* Can also be run as FIXED1

DRI Ethyl Glucuronide – Qualitative

Beckman Coulter System Parameters, AU2700/AU5400

Specific Test Parameters											
General		LIH	ISE	Range							
Test Name:	ETGQ		<	>	Type:	Urine		Operation:	Yes		
Sample:	Volume	25.0	μL	Dilution	0	μL	Pre-Dilution Rate:	1			
Reagents:	R1 Volume	57	μL	Dilution	0	μL	Min OD	Max OD			
	R2 Volume	57	μL	Dilution	0	μL	L	-2.00	H	3.00	
Wavelength:	Pri.	340	▽	Sec.	410	▽	Reagent OD limit:	First L	-2.00	First H	3.00
Method:	FIXED*				▽	Last L	-2.00	Last H	3.00		
Reaction slope:	+				▽	Dynamic Range:	L	#	H	#	
Measuring Point 1:	First	13	Last	17	Correlation Factor:	A	1	B	0		
Measuring Point 2:	First		Last		On-board stability period:	#					
Linearity					%						
No-Lag-Time:	No				▽						

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	ETGQ		<	>	Type:	Urine				
Value/Flag:	#		▽	Level L:	#		Level H:	#		
Normal Ranges:		Age L		Age H		L		H		
	Sex	Year	Month	Year	Month					
<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.	None Selected				#	#	#	#	#
	8.	Out of Range		L	H	#	#	#	#	#
Panic Value:	#		▽	#	Unit:	#		Decimal places:	#	

Calibration Specific										
General		ISE								
Test Name:	ETGQ		<	>	Type:	Urine				
Calibration Type:	AB		▽	Formula:	Y=AX+B		▽	Counts:	2	
	Process:	CONC								
	Cal. No.	OD	CONC	Factor/OD-L	Factor/OD-H					
Point 1:	#		100	-99999	99999					
Point 2:										
Point 3:										
Point 4:										
Point 5:										
Point 6:										
Point 7:										
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 FIXED

DRI Ethyl Glucuronide – Semiquantitative

Beckman Coulter System Parameters, AU2700/AU5400

Specific Test Parameters											
General		LIH	ISE	Range							
Test Name:	ETGSQ ▾			<	>	Type:	Urine ▾		Operation:	Yes ▾	
Sample:	Volume	25.0	μL	Dilution	0	μL	Pre-Dilution Rate:	1			
Reagents:	R1 Volume	57	μL	Dilution	0	μL	Min OD	Max OD			
	R2 Volume	57	μL	Dilution	0	μL	L	-2.00	H 2.50		
Wavelength:	Pri.	340	▾	Sec.	410	▾	Reagent OD limit:	First L	-2.00	First H	2.50
Method:	RATE1* ▾					Last L	-2.00	Last H	2.50		
Reaction slope:	+ ▾					Dynamic Range:	L	0	H 2000		
Measuring Point 1:	First	13		Last	17		Correlation Factor:	A	1	B	0
Measuring Point 2:	First			Last			On-board stability period:	#			
Linearity:						%					
No Lag Time:	No					▾					

Specific Test Parameters												
General		LIH	ISE	Range								
Test Name:	ETGSQ ▾			<	>	Type:	Urine ▾					
Value/Flag:	# ▾		Level L:	#		Level H:	#					
Normal Ranges:	Sex	Age L	Year	Month	Age H	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.	# ▾	#	#	#	#	#	#	#	#			
7. None Selected								#	#			
8. Out of Range								#	#			
Panic Value:			L			H			Unit:	ng/mL	Decimal places:	#

Calibration Specific										
General		ISE								
Test Name:	ETGSQ ▾			<	>	Type:	Urine ▾			
Calibration Type:	5AB ▾		Formula:	POLYGONAL ▾		Counts:	2		Process:	CONC ▾
Point 1:	Cal. No.	1	OD		CONC	0.0	Factor/OD-L	-2.0	Factor/OD-H	2.5
Point 2:	2			100.0	-2.0	2.5				
Point 3:	3			500.0	-2.0	2.5				
Point 4:	4			1000.0	-2.0	2.5				
Point 5:	5			2000.0	-2.0	2.5				
Point 6:										
Point 7:										
1-Point Cal. Point:				<input type="checkbox"/> With CONC-0						
MB Type Factor:				Calibration Stability Period:	#					

User defined
 * Can also be run as FIXED1

DRI Ethyl Glucuronide Assay – Qualitative

Beckman Coulter System Parameters, AU480 / AU680

Specific Test Parameters

General LIH ISE Range

Test Name: < > Type: Operation:

Sample Volume μ L Dilution μ L OD Limit

Pre-Dilution Rate Min. OD Max. OD

Reagents Volume: R1(R1-1) μ L Dilution μ L Reagent OD limit:

First Low High

Last Low High

R2 (R2-1) μ L Dilution μ L Dynamic Range Low High

Correlation Factor A B

Factor for Maker A B

Wavelength: Pri. nm Sec. nm

Method:

Reaction slope:

Onboard Stability Days Hour

Measuring Point 1: First Last LIH Influence Check

Measuring Point 2: First Last

Lipemia

Linearity: % Icterus

No Lag Time: Hemolysis

Specific Test Parameters

General ISE Range

Test Name: < > Type:

Value/Flag: Level L: Level H:

Specific Ranges:

	Sex	Year	Month	Year	Month	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="#"/>
<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="#"/>
<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="#"/>
<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="#"/>
<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="#"/>
<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="#"/>
7. No demographics						<input type="text" value="#"/>	<input type="text" value="#"/>
8. Not within expected values						<input type="text" value="#"/>	<input type="text" value="#"/>

Unit Decimal Places

Panic Value

Low High

Calibration Specific

General ISE

Test Name: < > Type: Use Serum Cal.

Calibration Type: Formula: Counts:

<Calibrator Parameters>

Calibrator †	OD	Conc	Factor Range		Slope Check <input type="text" value="+"/>
			Low	High	
Point 1: <input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="100"/>	<input type="text" value="-99999"/>	<input type="text" value="999999"/>	Allowable Range Check <input type="checkbox"/> Reagent Blank <input type="checkbox"/> Calibration Advanced Calibration Operation <input type="text" value="No"/> Interval (RB/ACAL) <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=""/>	
Point 3: <input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point 4: <input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point 5: <input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point 6: <input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point 7: <input type="text" value=""/>	<input type="text" value=""/>	<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=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point 9: <input type="text" value=""/>	<input type="text" value=""/>	<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=""/>	<input type="text" value=""/>	<input type="text" value=""/>	

<Point Cal. For Master Curve>

Calibrator	OD	Conc	OD Range		Stability
			Low	High	
Point 1: <input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Reagent Blanks Calibration <input type="text" value="#"/> Day <input type="text" value="30"/> Hour
Point 2: <input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Reagent Blanks Calibration <input type="text" value="#"/> Day <input type="text" value="0"/> Hour

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

Lot Calibration

User defined.

*Can also be run as RATE

DRI Ethyl Glucuronide Assay – Semiquantitative Beckman Coulter System Parameters, AU480 /AU680

Specific Test Parameters

General | LIH | ISE | Range

Test Name: < > Type: Operation:

Sample Volume μL Dilution μL OD Limit

Pre-Dilution Rate Min. OD Max. OD

Reagents Volume: R1(R1-1) μL Dilution μL Reagent OD limit:

First Low Hig

Last Low Hig

R2 (R2-1) μL Dilution μL Dynamic Range Low Hig

Correlation Factor A B

Factor for Maker A B

Wavelength: Pri. nm Sec. nm

Method:

Reaction slope:

Onboard Stability Days Hour

Measuring Point 1: First Last LIH Influence Check

Measuring Point 2: First Last

Lipemia

Linearity: % Icterus

No Lag Time: Hemolysis

Specific Test Parameters

General | ISE | Range

Test Name: < > Type:

Value/Flag: Level L: Level H:

Specific Ranges:

	Sex	Year	Month	Year	Month	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="#"/>
<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="#"/>
<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="#"/>
<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="#"/>
<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="#"/>
<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="#"/>
7. No demographics						<input type="text" value="#"/>	<input type="text" value="#"/>
8. Not within expected values						<input type="text" value="#"/>	<input type="text" value="#"/>

Unit Decimal Places

Panic Value

Low High

Calibration Specific

General | ISE

Test Name: < > Type: Use Serum Cal.

Calibration Type: Formula: Counts:

<Calibrator Parameters> Slope Check

Calibrator †	OD	Conc	Factor Range		Allowable Range Check
			Low	High	
Point 1:	<input type="text" value="#"/>	<input type="text" value="0.0"/>	<input type="text" value="-2.00"/>	<input type="text" value="3.00"/>	<input type="checkbox"/> Reagent Blank
Point 2:	<input type="text" value="#"/>	<input type="text" value="100.0"/>	<input type="text" value="-2.00"/>	<input type="text" value="3.00"/>	<input type="checkbox"/> Calibration
Point 3:	<input type="text" value="#"/>	<input type="text" value="500.0"/>	<input type="text" value="-2.00"/>	<input type="text" value="3.00"/>	Advanced Calibration
Point 4:	<input type="text" value="#"/>	<input type="text" value="1000.0"/>	<input type="text" value="-2.00"/>	<input type="text" value="3.00"/>	
Point 5:	<input type="text" value="#"/>	<input type="text" value="2000.0"/>	<input type="text" value="-2.00"/>	<input type="text" value="3.00"/>	Operation <input type="text" value="No"/>
Point 6:	<input type="text" value=""/>				Interval (RB/ACAL) <input type="text" value=""/>
Point 7:	<input type="text" value=""/>				
Point 8:	<input type="text" value=""/>				
Point 9:	<input type="text" value=""/>				
Point 10:	<input type="text" value=""/>				

<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:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value="#"/> Day <input type="text" value="#"/> Hour
Point 2:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value="30"/> Day <input type="text" value="0"/> Hour

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

User Defined
* Can also be run as RATE1

DRI Ethyl Glucuronide Assay – Qualitative

Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: <input type="text" value="ETGQ"/> < > Type: <input type="text" value="Urine"/> Operation <input type="text" value="Yes"/>					
Sample Volume	<input type="text" value="17.0"/> μL	Dilution	<input type="text" value="0"/> μL	OD Limit	
Pre-Dilution Rate	<input type="text" value="1"/>	Diluent Bottle	<input type="text" value="Outside"/>	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="39"/> μL	Dilution	<input type="text" value="0"/> μL	Reagent OD Limit	
	R1-2 <input type="text" value=""/>	Dilution	<input type="text" value=""/>	1 st . Low <input type="text" value="-2.00"/> High <input type="text" value="3.00"/>	
	R2(R2-1) <input type="text" value="39"/> μL	Dilution	<input type="text" value="0"/> μL	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="None"/>	Dynamic Range Low	<input type="text" value="#"/>
Wavelength	Pri <input type="text" value="340"/> nm	Sec.	<input type="text" value="410"/> nm	Correlation Factor A	<input type="text" value="1"/>
Method	<input type="text" value="FIXED*"/>			Factor for Maker A	<input type="text" value="1"/>
Reaction Slope	<input type="text" value="+"/>			High B	<input type="text" value="0"/>
Measuring Point1 1 st	<input type="text" value="13"/>	Last	<input type="text" value="17"/>	Onboard Stability Period	<input type="text" value="30"/> Day <input type="text" value="0"/> Hour
Measuring Point2 1 st	<input type="text" value=""/>	Last	<input type="text" value=""/>	LIH Influence Check	<input type="text" value="No"/>
Linearity Limit	<input type="text" value=""/>			Lipemia	<input type="text" value=""/>
Lag Time Check	<input type="text" value="No"/>			Icterus	<input type="text" value=""/>
				Hemolysis	<input type="text" value=""/>

Parameters		Specific Test Parameters						
General	LIH	ISE	HbA1c	Calculated Test	Range			
Test Name: <input type="text" value="ETGQ"/> < > Type: <input type="text" value="Urine"/>								
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	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	<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="ETGQ"/> < > Type: <input type="text" value="Urine"/> Cuvette: <input type="text" value=""/>					
<input type="checkbox"/> Use Serum Cal.					
Calibration Type: <input type="text" value="AB"/> Formula: <input type="text" value="Y=AX+B"/> Counts: <input type="text" value="2"/>					
<Calibrator Parameters> Range					
Calibrator	OD	Conc	Low	High	Slope Check
Point 1:	<input type="text" value="#"/>	<input type="text" value="100"/>	<input type="text" value="-99999"/>	<input type="text" value="999999"/>	<input type="text" value="+"/>
Point 2:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Allowance Range Check
Point 3:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="checkbox"/> Reagent Blank
Point 4:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="checkbox"/> Calibration
Point 5:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Advanced Calibration
Point 6:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Operation
Point 7:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value="No"/>
Point 8:	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Interval (RB/ACAL)
Point 9:	<input type="text" value=""/>	<input type="text" value=""/>	<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=""/>	<input type="text" value=""/>	<input type="checkbox"/> Lot Calibration
<Point Cal. For No. of Correction Points <input type="text" value=""/> Use Master Curve <input type="text" value=""/>					
Master Curve> OD Range					
Calibrator	OD	Conc	Low	High	Stability
Point-1	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Reagent Blank
Point-2	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Calibration
MB Type Factor: <input type="text" value=""/> 1-Point Calibration Point <input type="text" value="None"/> <input type="checkbox"/> with Conc-0					

User defined.
* Can also be run as RATE

DRI Ethyl Glucuronide Assay – Semiquantitative Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters					
General		LIH	ISE	HbA1c		Calculated Test	Range
Test Name:		ETGSQ		<	>	Type:	Urine
						Operation	Yes
Sample Volume	17.0	μL	Dilution	0	μL	OD Limit	
Pre-Dilution Rate	1	▽	Diluent Bottle	Outside		Min.OD	-2.00
Rgt. Volume	39	μL	Dilution	0	μL	Max.OD	3.00
	R1(R1-1)		Dilution			Reagent OD Limit	
	R1-2		Dilution			1 st	Low -2.00 High 3.00
	R2(R2-1)	39	μL	Dilution	0	Last	Low -2.00 High 3.00
Common Rgt. Type	None		Name	None		Dynamic Range Low	0
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 Period	30 Day 0 Hour
Measuring Point1 1 st	13		Last	17		LIH Influence Check	No
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:		ETGSQ		<	>	Type:	Urine	
Value/Flag:		#						
						Level	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.	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:		ETGSQ		<	>	Type:	Urine
						Cuvette .	
						<input type="checkbox"/> Use Serum Cal.	
Calibration Type:		5AB		Formula:	Polygonal		Counts: 2
<Calibrator Parameters>				Range			
Calibrator	OD	Conc	Low	High	Slope Check		
Point 1:	#	0.0	-2.00	3.00	+		
Point 2:	#	100.0	-2.00	3.00	Allowance Range Check		
Point 3:	#	500.0	-2.00	3.00	<input type="checkbox"/> Reagent Blank		
Point 4:	#	1000.0	-2.00	3.00	<input type="checkbox"/> Calibration		
Point 5:	#	2000.0	-2.00	3.00	Advanced Calibration		
Point 6:					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>							
Calibrator	OD	Conc	Low	High	Stability		
Point-1					Reagent Blank # Day # Hour		
Point-2					Calibration 30 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:

Qualitative Assay 500 ng/mL Cutoff

Controls	Low Control	Cutoff Cal	High Control
AU480 (N=84)			
Mean Rate (mA/min)	436.8	457.1	480.1
Within-Run SD (mA/min)	1.96	1.35	2.12
Within-Run CV (%)	0.5	0.3	0.4
Total SD (mA/min)	2.64	2.06	2.73
Total CV (%)	0.6	0.5	0.6
AU640 (N=60)			
Mean Rate (mA/min)	382	401	423
Within-Run SD (mA/min)	2.6	1.6	3.3
Within-Run CV (%)	0.7	0.4	0.8
Total SD (mA/min)	4.1	5.5	6.0
Total CV (%)	1.1	1.4	1.4
AU680 (N=84)			
Mean Rate (mA/min)	455.4	476.1	503.4
Within-Run SD (mA/min)	1.14	1.40	1.35
Within-Run CV (%)	0.3	0.3	0.3
Total SD (mA/min)	2.59	3.10	3.56
Total CV (%)	0.6	0.7	0.7
AU5800 (N=84)			
Mean Rate (mA/min)	438.7	457.0	480.4
Within-Run SD (mA/min)	1.40	1.46	2.01
Within-Run CV (%)	0.3	0.3	0.4
Total SD (mA/min)	2.54	2.87	3.27
Total CV (%)	0.6	0.6	0.7

**Precision,
continued**

**Qualitative Assay
1000 ng/mL Cutoff**

Controls	Low Control	Cutoff Cal	High Control
AU480 (N=80)			
Mean Rate (mA/min)	530.0	561.2	590.8
Within-Run SD (mA/min)	2.64	2.58	2.53
Within-Run CV (%)	0.5	0.5	0.4
Total SD (mA/min)	2.85	2.78	2.83
Total CV (%)	0.5	0.5	0.5
AU640 (N=60)			
Mean Rate (mA/min)	450	474	506
Within-Run SD (mA/min)	4.6	3.2	4.7
Within-Run CV (%)	1.0	2.2	0.9
Total SD (mA/min)	6.0	10.5	7.1
Total CV (%)	1.3	2.2	1.4
AU680 (N=80)			
Mean Rate (mA/min)	557.6	589.3	620.4
Within-Run SD (mA/min)	1.90	3.00	2.93
Within-Run CV (%)	0.3	0.5	0.5
Total SD (mA/min)	2.96	4.52	4.07
Total CV (%)	0.5	0.8	0.7
AU5800 (N=84)			
Mean Rate (mA/min)	524.4	554.2	584.2
Within-Run SD (mA/min)	4.95	5.05	4.91
Within-Run CV (%)	0.9	0.9	0.8
Total SD (mA/min)	7.26	7.30	7.59
Total CV (%)	1.4	1.3	1.3

**Precision,
continued**

**Semiquantitative Assay
500 ng/mL Cutoff**

Controls	Low Control	Cutoff Cal	High Control
AU480 (N=84)			
Mean (ng/mL)	407.6	507.1	646.4
Within-Run SD (ng/mL)	9.52	7.94	12.91
Within-Run CV (%)	2.3	1.6	2.0
Total SD (ng/mL)	11.99	8.76	13.31
Total CV (%)	2.9	1.7	2.1
AU640 (N=60)			
Mean (ng/mL)	382	486	612
Within-Run SD (ng/mL)	17.0	9.7	21.6
Within-Run CV (%)	4.5	2.0	3.5
Total SD (ng/mL)	21.1	25.9	30.9
Total CV (%)	5.5	5.3	5.1
AU680 (N=84)			
Mean (ng/mL)	400.8	489.6	647.1
Within-Run SD (ng/mL)	4.89	6.37	8.04
Within-Run CV (%)	1.2	1.3	1.2
Total SD (ng/mL)	9.47	10.89	12.08
Total CV (%)	2.4	2.2	1.9
AU5800 (N=84)			
Mean (ng/mL)	412.3	499.7	639.5
Within-Run SD (ng/mL)	6.65	7.71	12.13
Within-Run CV (%)	1.6	1.5	1.9
Total SD (ng/mL)	7.88	11.20	14.58
Total CV (%)	1.9	2.2	2.3

**Precision,
continued**

**Semiquantitative Assay
1000 ng/mL Cutoff**

Controls	Low Control	Cutoff Cal	High Control
AU480 (N=80)			
Mean (ng/mL)	777.1	1007.8	1244.3
Within-Run SD (ng/mL)	19.40	19.96	20.17
Within-Run CV (%)	2.5	2.0	1.6
Total SD (ng/mL)	20.15	23.13	22.76
Total CV (%)	2.6	2.3	1.8
AU640 (N=60)			
Mean (ng/mL)	810	996	1291
Within-Run SD (ng/mL)	26.8	25.8	39.5
Within-Run CV (%)	3.3	2.6	3.1
Total SD (ng/mL)	37.6	57.6	59.2
Total CV (%)	4.7	5.8	4.6
AU680 (N=80)			
Mean (ng/mL)	781.3	1004.6	1241.7
Within-Run SD (ng/mL)	13.42	21.63	21.92
Within-Run CV (%)	1.7	2.2	1.8
Total SD (ng/mL)	21.15	33.19	31.78
Total CV (%)	2.7	3.3	2.6
AU5800 (N=84)			
Mean (ng/mL)	771.3	990.4	1236.7
Within-Run SD (ng/mL)	28.99	21.56	31.23
Within-Run CV (%)	3.8	2.2	2.5
Total SD (ng/mL)	38.01	39.00	44.46
Total CV (%)	4.9	3.9	3.6

Linearity

Nine levels of diluted calibrators were run against a single calibration curve and the linearity calculated.

The Ethyl Glucuronide assay recovered between 99.0% and 106.4% on the AU480.
The Ethyl Glucuronide assay recovered between 93.2% and 106.3% on the AU680.
The Ethyl Glucuronide assay recovered between 98.5% and 108.3% on the AU5800.

**Limit of
Detection
(LDD)**

The negative calibrator was run against the same calibration curve for 21 replicates.
The LDD is calculated as 2*SD.

The observed LDDs for DRI Ethyl Glucuronide on the AU480, AU680 and AU5800 were 13.5, 9.7, and 11.4 ng/mL, respectively.

Accuracy and Correlation

Qualitative 500 ng/mL

Patient urine samples were assayed using the DRI Ethyl Glucuronide Assay on the Beckman Coulter AU480, AU640, AU680, and AU5800 and tested against the reference analyzer, the Hitachi 717.

Qualitative (500 ng/mL Cutoff) Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	94.3%	100.0%	96.2%
AU640	97.9%	100.0%	99.0%
AU680	94.3%	100.0%	96.2%
AU5800	97.7%	100.0%	98.5%

DRI Ethyl Glucuronide Assay Qualitative (500 ng/mL) Method Comparison

		Hitachi 717	
		+	-
AU480	+	83	0
	-	5	42

		Hitachi 717	
		+	-
AU640	+	46	0
	-	1	53

		Hitachi 717	
		+	-
AU680	+	83	0
	-	5	42

		Hitachi 717	
		+	-
AU5800	+	86	0
	-	2	42

Accuracy and Correlation

Qualitative 1000 ng/mL

Patient urine samples were assayed using the DRI Ethyl Glucuronide Assay on the Beckman Coulter AU480, AU640, AU680, and AU5800 and tested against the reference analyzer, the Hitachi 717.

Qualitative (1000 ng/mL Cutoff) Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	98.5%	100.0%	99.2%
AU640	93.8%	100.0%	98.0%
AU680	95.4%	100.0%	97.7%
AU5800	98.5%	96.9%	97.7%

DRI Ethyl Glucuronide Assay Qualitative (1000 ng/mL) Method Comparison

		Hitachi 717	
		+	-
AU480	+	64	0
	-	1	65

		Hitachi 717	
		+	-
AU640	+	30	0
	-	2	68

		Hitachi 717	
		+	-
AU680	+	62	0
	-	3	65

		Hitachi 717	
		+	-
AU5800	+	64	2
	-	1	63

Accuracy and Correlation

Semiquantitative 500 ng/mL

Patient urine samples were assayed using the DRI Ethyl Glucuronide Assay on the Beckman Coulter AU480, AU640, AU680, and AU5800 and tested against the reference analyzer, the Hitachi 717.

Semiquantitative (500 ng/mL Cutoff) Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	98.8%	100.0%	99.2%
AU640	100.0%	100.0%	100.0%
AU680	100.0%	97.9%	99.2%
AU5800	100.0%	100.0%	100.0%

**DRI Ethyl Glucuronide Assay
Semiquantitative (500 ng/mL) Method Comparison**

		Hitachi 717				Hitachi 717	
		+	-			+	-
AU480	+	83	0	AU640	+	46	0
	-	1	46		-	0	54
		Hitachi 717				Hitachi 717	
		+	-			+	-
AU680	+	83	1	AU5800	+	84	0
	-	0	46		-	0	46

Accuracy and Correlation

Semiquantitative 1000 ng/mL

Patient urine samples were assayed using the DRI Ethyl Glucuronide Assay on the Beckman Coulter AU480, AU640, AU680, and AU5800 and tested against the reference analyzer, the Hitachi 717.

Semiquantitative (1000 ng/mL Cutoff) Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	98.4%	98.5%	98.5%
AU680	100.0%	98.5%	99.2%
AU5800	98.4%	100.0%	99.2%

**DRI Ethyl Glucuronide Assay
Semiquantitative (1000 ng/mL) Method Comparison**

		Hitachi 717	
		+	-
AU480	+	63	1
	-	1	65

		Hitachi 717	
		+	-
AU680	+	63	1
	-	0	66

		Hitachi 717	
		+	-
AU5800	+	63	0
	-	1	66

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.

End