

DRI[®] ETHYL GLUCURONIDE APPLICATION BECKMAN COULTER DxC 700 AU[®]



Beckman Coulter Reagent REF E0750310

Intended for the qualitative and semi-quantitative determination of Ethyl Glucuronide in human urine at cutoffs of 500 ng/mL and 1000 ng/mL. This application pertains only to the 500 ng/mL cutoff.



For In Vitro Diagnostic Use Only
For Export Only, Not Available in the United States
Rx 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
DRI Ethyl Glucuronide Assay	1 x 18 mL	E0750310
DRI Ethyl Glucuronide Negative Calibrator	1 x 25 mL	E0750311
DRI Ethyl Glucuronide 100 Calibrator	1 x 10 mL	E0750312
DRI Ethyl Glucuronide 500 Calibrator	1 x 10 mL	E0750313
DRI Ethyl Glucuronide 1000 Calibrator	1 x 10 mL	E0750314
DRI Ethyl Glucuronide 2000 Calibrator	1 x 10 mL	E0750315
DRI Ethyl Glucuronide 375 Control	1 x 25 mL	E0750316
DRI Ethyl Glucuronide 625 Control	1 x 25 mL	E0750317
AU Bottle	20 X 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. Refer to the package insert for complete reagent preparation.

Prior to pouring into AU bottles, allow the reagent to equilibrate for 15 minutes at refrigerated temperature (2 to 8°C). Dispense R1 reagent and R2 reagent into appropriate AU bottles as shown in the table below:

DRI Ethyl Glucuronide Assay Kit	AU Reagent Bottle	
	R1 Compartment	R2 Compartment
Antibody/Substrate Reagent <u>R1</u>	One Bottle (30 mL)	
Enzyme Conjugate Reagent <u>R2</u>		One Bottle (30 mL)

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

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. The product insert can be found at the Thermo Fisher Scientific website:

www.thermofisher.com

Calibration

Use the DRI Ethyl Glucuronide calibrators. The calibrators are liquid and ready-to-use.



DRI Ethyl Glucuronide Assay
Beckman Coulter System Parameters, DxC 700 AU - Qualitative

General	LIH	ISE	Calculated Test	Range
Test Name: # <input type="text"/> Type: Urine <input type="text"/> Operation: Yes <input type="text"/>				
Sample Volume	<input type="text" value="25.0"/> μL	Dilution	<input type="text" value="0"/> μL	OD Limit
Pre-Dilution Rate	<input type="text" value="1"/>			Min. OD <input type="text" value="-2.0000"/> Max OD <input type="text" value="3.0000"/>
Reagent Volume	R1 (R1-1) <input type="text" value="57"/> μL	Dilution	<input type="text" value="0"/> μL	Reagent OD Limit 1 st Low <input type="text" value="-2.0000"/> High <input type="text" value="3.0000"/>
	R1-2 <input type="text"/> μL	Dilution	<input type="text"/> μL	Last Low <input type="text" value="-2.0000"/> High <input type="text" value="3.0000"/>
	R2 (R2-1) <input type="text" value="57"/> μL	Dilution	<input type="text" value="0"/> μL	Analytical Measuring Range Low <input type="text" value="-9999"/> High <input type="text" value="9999"/>
Common Reagent	Type <input type="text" value="None"/>	Name	<input type="text" value="None"/>	Correlation Factor A <input type="text" value="1"/> B <input type="text" value="0"/>
Wavelength	<input type="text" value="340"/> nm	Sec	<input type="text" value="410"/> nm	Manufacturer Factor A <input type="text" value="1"/> B <input type="text" value="0"/>
Method	<input type="text" value="FIXED"/>			Onboard Stability Period <input type="text" value="30"/> Day <input type="text" value="0"/> Hour
Reaction Slope	<input type="text" value="+"/> \downarrow			LIH Influence Check <input type="text" value="No"/> \downarrow
Measuring Point-1	1st <input type="text" value="13"/>	Last	<input type="text" value="17"/>	Lipemia <input type="text" value="+"/> \downarrow
Measuring Point-2	1st <input type="text"/>	Last	<input type="text"/>	Icterus <input type="text" value="+"/> \downarrow
Linearity Limit	<input type="text"/> %			Hemolysis <input type="text" value="+"/> \downarrow
Lag Time Check	<input type="text" value="No"/> \downarrow			

General	LIH	ISE	Calculated Test	Range
Test Name: # <input type="text"/> Type: Urine <input type="text"/>				
Value/Flag	<input type="text" value="Flag"/>	Level	Low <input type="text" value="-999999"/> High <input type="text" value="500"/>	
Specific Ranges				
	Sex	From	To	Other Type
	Year	Month	Year	Month
<input type="checkbox"/> 1:	<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="checkbox"/> 3:	<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="checkbox"/> 5:	<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="#"/>
7:	Standard demographics			<input type="text" value="#"/>
8:	Not within expected values			<input type="text" value="#"/>
Critical Limits	Low <input type="text" value="#"/>	High <input type="text" value="#"/>	Unit <input type="text" value="ng/mL"/>	Select <input type="text" value="0"/> Decimal Places

User-defined



DRI Ethyl Glucuronide Assay
Beckman Coulter System Parameters, DxC 700 AU – Qualitative, *continued*

Calibrators	General	ISE																																																			
Test Name: # Type: Urine																																																					
<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> <table border="1" style="width:100%; border-collapse: collapse; margin-top: 10px;"> <thead> <tr> <th rowspan="2"></th> <th rowspan="2">Calibrator</th> <th rowspan="2">OD</th> <th rowspan="2">Conc</th> <th colspan="2">Range</th> </tr> <tr> <th>Low</th> <th>High</th> </tr> </thead> <tbody> <tr> <td>Point-1</td> <td>#</td> <td></td> <td>500</td> <td>-999999</td> <td>999999</td> </tr> <tr> <td>Point-2</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> <tr> <td>Point-3</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> <tr> <td>Point-4</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> <tr> <td>Point-5</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> <tr> <td>Point-6</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> <tr> <td>Point-7</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>					Calibrator	OD	Conc	Range		Low	High	Point-1	#		500	-999999	999999	Point-2						Point-3						Point-4						Point-5						Point-6						Point-7					
	Calibrator	OD	Conc					Range																																													
				Low	High																																																
Point-1	#		500	-999999	999999																																																
Point-2																																																					
Point-3																																																					
Point-4																																																					
Point-5																																																					
Point-6																																																					
Point-7																																																					
MB Type Factor <input type="text"/>		1-Point Calibration Point <input type="text" value="None"/>																																																			
		<input type="checkbox"/> with Conc-0 <table style="margin-left: 20px;"> <tr> <td>Slope Check</td> <td><input type="text" value="1"/></td> </tr> <tr> <td>Allowable Range Check</td> <td><input type="text"/></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)</td> <td><input type="text"/></td> </tr> <tr> <td>Interval (ACAL)</td> <td><input type="text"/></td> </tr> <tr> <td>Stability</td> <td></td> </tr> <tr> <td>Reagent Blank</td> <td><input type="text" value="#"/> Day <input type="text" value="0"/> Hour</td> </tr> <tr> <td>Calibration</td> <td><input type="text" value="#"/> Day <input type="text" value="0"/> Hour</td> </tr> </table>		Slope Check	<input type="text" value="1"/>	Allowable Range Check	<input type="text"/>	<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)	<input type="text"/>	Interval (ACAL)	<input type="text"/>	Stability		Reagent Blank	<input type="text" value="#"/> Day <input type="text" value="0"/> Hour	Calibration	<input type="text" value="#"/> Day <input type="text" value="0"/> Hour																														
Slope Check	<input type="text" value="1"/>																																																				
Allowable Range Check	<input type="text"/>																																																				
<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)	<input type="text"/>																																																				
Interval (ACAL)	<input type="text"/>																																																				
Stability																																																					
Reagent Blank	<input type="text" value="#"/> Day <input type="text" value="0"/> Hour																																																				
Calibration	<input type="text" value="#"/> Day <input type="text" value="0"/> Hour																																																				

User-defined



DRI Ethyl Glucuronide Assay
Beckman Coulter System Parameters, DxC 700 AU – Semiquantitative

General	LIH	ISE	Calculated Test	Range
Test Name: # <input type="text"/> Type: Urine Operation: Yes				
Sample Volume	<input type="text" value="25.0"/> μL	Dilution <input type="text" value="0"/> μL	OD Limit	
Pre-Dilution Rate	<input type="text" value="1"/>		Min. OD <input type="text" value="-2.0000"/> Max OD <input type="text" value="3.0000"/>	
Reagent Volume	R1 (R1-1) <input type="text" value="57"/> μL	Dilution <input type="text" value="0"/> μL	Reagent OD Limit 1 st	Low <input type="text" value="-2.0000"/> High <input type="text" value="3.0000"/>
	R1-2 <input type="text"/> μL	Dilution <input type="text"/> μL	Last	Low <input type="text" value="-2.0000"/> High <input type="text" value="3.0000"/>
	R2 (R2-1) <input type="text" value="57"/> μL	Dilution <input type="text" value="0"/> μL	Analytical Measuring Range	Low <input type="text" value="-9999"/> High <input type="text" value="9999"/>
Common Reagent	Type <input type="text" value="None"/>	Name <input type="text" value="None"/>	Correlation Factor	A <input type="text" value="1"/> B <input type="text" value="0"/>
Wavelength	<input type="text" value="340"/> nm	Sec <input type="text" value="410"/> nm	Manufacturer Factor	A <input type="text" value="1"/> B <input type="text" value="0"/>
Method	<input type="text" value="FIXED1"/>		Onboard Stability Period	<input type="text" value="30"/> Day <input type="text" value="0"/> Hour
Reaction Slope	<input type="text" value="+"/> μV		LIH Influence Check	<input type="text" value="No"/>
Measuring Point-1	1st <input type="text" value="13"/> Last <input type="text" value="17"/>		Lipemia	<input type="text" value="+"/> μV
Measuring Point-2	1st <input type="text"/> Last <input type="text"/>		Icterus	<input type="text" value="+"/> μV
Linearity Limit	<input type="text"/> %		Hemolysis	<input type="text" value="+"/> μV
Lag Time Check	<input type="text" value="No"/>			

General	LIH	ISE	Calculated Test	Range
Test Name: # <input type="text"/> Type: Urine				
Value/Flag	<input type="text" value="Value"/>	Level	Low <input type="text" value="-999999"/> High <input type="text" value="500"/>	
Specific Ranges				
	Sex	From Year	To Year	Other Type
<input type="checkbox"/> 1:	<input type="text" value="#"/>	Month <input type="text" value="#"/>	Month <input type="text" value="#"/>	None <input type="text" value="None"/>
<input type="checkbox"/> 2:	<input type="text" value="#"/>	Month <input type="text" value="#"/>	Month <input type="text" value="#"/>	None <input type="text" value="None"/>
<input type="checkbox"/> 3:	<input type="text" value="#"/>	Month <input type="text" value="#"/>	Month <input type="text" value="#"/>	None <input type="text" value="None"/>
<input type="checkbox"/> 4:	<input type="text" value="#"/>	Month <input type="text" value="#"/>	Month <input type="text" value="#"/>	None <input type="text" value="None"/>
<input type="checkbox"/> 5:	<input type="text" value="#"/>	Month <input type="text" value="#"/>	Month <input type="text" value="#"/>	None <input type="text" value="None"/>
<input type="checkbox"/> 6:	<input type="text" value="#"/>	Month <input type="text" value="#"/>	Month <input type="text" value="#"/>	None <input type="text" value="None"/>
7:	Standard demographics			Low <input type="text" value="#"/> High <input type="text" value="#"/>
8:	Not within expected values			Low <input type="text" value="#"/> High <input type="text" value="#"/>
Critical Limits	Low <input type="text" value="#"/>	High <input type="text" value="#"/>	Unit <input type="text" value="ng/mL"/>	Decimal Places <input type="text" value="1"/>

User-defined



**DRI Ethyl Glucuronide Assay
Beckman Coulter System Parameters, DxC 700 AU – Semiquantitative, *continued***

Calibrators	General	ISE																																																		
Test Name: # Type: Serum																																																				
<input type="checkbox"/> Use Serum Cal.																																																				
Calibration Type: 5AB Formula: POLYGONAL Counts: 2																																																				
<Calibrator Parameters>																																																				
<table border="1"> <thead> <tr> <th rowspan="2"></th> <th rowspan="2">Calibrator</th> <th rowspan="2">OD</th> <th rowspan="2">Conc</th> <th colspan="2">Range</th> </tr> <tr> <th>Low</th> <th>High</th> </tr> </thead> <tbody> <tr> <td>Point-1</td> <td>#</td> <td></td> <td>0</td> <td>-2.00</td> <td>3.00</td> </tr> <tr> <td>Point-2</td> <td>#</td> <td></td> <td>100</td> <td>-2.00</td> <td>3.00</td> </tr> <tr> <td>Point-3</td> <td>#</td> <td></td> <td>500</td> <td>-2.00</td> <td>3.00</td> </tr> <tr> <td>Point-4</td> <td>#</td> <td></td> <td>1000</td> <td>-2.00</td> <td>3.00</td> </tr> <tr> <td>Point-5</td> <td>#</td> <td></td> <td>2000</td> <td>-2.00</td> <td>3.00</td> </tr> <tr> <td>Point-6</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> <tr> <td>Point-7</td> <td></td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>				Calibrator	OD	Conc	Range		Low	High	Point-1	#		0	-2.00	3.00	Point-2	#		100	-2.00	3.00	Point-3	#		500	-2.00	3.00	Point-4	#		1000	-2.00	3.00	Point-5	#		2000	-2.00	3.00	Point-6						Point-7					
	Calibrator	OD					Conc	Range																																												
			Low	High																																																
Point-1	#		0	-2.00	3.00																																															
Point-2	#		100	-2.00	3.00																																															
Point-3	#		500	-2.00	3.00																																															
Point-4	#		1000	-2.00	3.00																																															
Point-5	#		2000	-2.00	3.00																																															
Point-6																																																				
Point-7																																																				
MB Type Factor: 1-Point Calibration Point: None																																																				
<input type="checkbox"/> with Conc=0																																																				
Slope Check: +																																																				
<input type="checkbox"/> Allowable Range Check																																																				
<input type="checkbox"/> Reagent Blank																																																				
<input type="checkbox"/> Calibration																																																				
Advanced Calibration Operation: No																																																				
Interval (RB):																																																				
Interval (ACAL):																																																				
Stability																																																				
Reagent Blank: # Day 0 Hour																																																				
Calibration: # Day 0 Hour																																																				

User-defined

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 Technical Support Center if this product is received damaged.

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AU Series Systems are trademarks of Beckman Coulter.



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

End