

DRI™ ETHYL ALCOHOL ASSAY APPLICATION CE
BECKMAN COULTER AU400® , AU480® , AU640® , AU680® ,
AU2700® , AU5400® , AU5800®

Catalog No. 0037, 0038, 10016397

This Application is Intended for the Quantitative Determination of Alcohol in Human Urine, Serum or Plasma



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 manual for information on analyzer operation.

Dispense adequate amounts of Buffer Reagent A (first reagent) and Enzyme 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.**

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

Quantitative results will be printed in mg/dL.

DRI Ethyl Alcohol Assay - Quantitative

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

Specific Test Parameters	
General	LIH ISE Range
Test Name:	# < > Type: # Operation: Yes
Sample:	Volume: 9.0 μ L Dilution: 0 μ L Pre-Dilution Rate: 1
Reagents:	R1 Volume: 75 μ L Dilution: 0 μ L Min OD: Max OD
	R2 Volume: 75 μ L Dilution: 0 μ L L: -2.00 H: 2.50
Wavelength:	Pri.: 340 μ L Sec.: 410 μ L Reagent OD limit: First L: -2.00 First H: 2.50
Method:	RATE1* Last L: -2.00 Last H: 2.50
Reaction slope:	+
Measuring Point 1:	First: 13 Last: 17 Dynamic Range: L: 10 H: 600
Measuring Point 2:	First: Last: Correlation Factor: A: 1 B: 0
Linearity:	% On-board stability period: #
No Lag Time:	No

General		LIH	ISE	Range	
Test Name:	# < >	Type:	#		
Value/Flag:	#	Level L:	#	Level H:	#
Normal Ranges:	Sex	Age L	Age H	L	H
	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				-99999	100
8. Out of Range		L	H	#	#
Panic Value:	#	#	#	Unit: mg/dL	Decimal places: #

Calibration Specific	
General	ISE
Test Name:	# < > Type: #
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: 999999
Point 2:	Cal. No.: # OD: CONC: 100£ Factor/OD-L: -99999 Factor/OD-H: 999999
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
MB Type Factor:	Calibration Stability Period: #

User Defined

* Can be run as FIXED1

£ For conversion to g/L change the calibrator concentration to 1 and the units to g/L

NOTE : If your software has a slope check field, then input "+" into that field

DRI Ethyl Alcohol Assay - Quantitative

Beckman Coulter System Parameters, AU480/AU680

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	#	<	>	Type:	#	Operation:	Yes			
Sample Volume	9.0	μL	Dilution	0	μL	OD Limit				
Pre-Dilution Rate	1					Min. OD	-2.00	Max. OD	3.00	
Reagents Volume:	R1(R1-1)	75	μL	Dilution	0	μL	Reagent OD limit:			
						First Low	-2.00	Hig	3.00	
						Last Low	-2.00	Hig	3.00	
	R2 (R2-1)	75	μL	Dilution	0	μL	Dynamic Range Low	10	Hig	600
						Correlation Factor A	1	B	0	
Wavelength:	Pri.	340	nm	Sec.	410	nm	Factor for Maker A	1	B	0
Method:	FIXED1* ▾									
Reaction slope:	+ ▾									
Measuring Point 1:	First	13	Last	17	Onboard Stability <input type="checkbox"/> Days <input type="checkbox"/> Hour					
Measuring Point 2:	First		Last		LIH Influence Check # ▾					
Linearity:	% ▾									
No Lag Time:	No ▾									
						Lipemia	▾			
						Icterus	▾			
						Hemolysis	▾			

Specific Test Parameters										
General		ISE	Range							
Test Name:	#	<	>	Type:	#					
Value/Flag:	VALUE	Level L:	#	Level H:	#	Panic Value				
Specific Ranges:		From	To		Low	High				
<input type="checkbox"/>	1.	Sex	Year	Month	Year	Month	Low	High	Low High	
<input type="checkbox"/>	2.	#	#	#	#	#	#	#	# #	
<input type="checkbox"/>	3.	#	#	#	#	#	#	#	# #	
<input type="checkbox"/>	4.	#	#	#	#	#	#	#	# #	
<input type="checkbox"/>	5.	#	#	#	#	#	#	#	# #	
<input type="checkbox"/>	6.	#	#	#	#	#	#	#	# #	
<input type="checkbox"/>	7.	No demographics					#	#		
<input type="checkbox"/>	8.	Not within expected values					#	#		
	Unit	mg/dL		Decimal Places	#					

* Can be run as RATE1
User Defined Values

DRI Ethyl Alcohol Assay - Quantitative
Beckman Coulter System Parameters, AU480/AU680, *continued*

Calibration Specific									
General		ISE							
Test Name:		#	<	>	Type	#	<input type="checkbox"/> Use Serum Cal.		
Calibration Type:		AA	Formula:		Y=Ax+B	Counts:	2		
<Calibrator Parameters>									
	Calibrator †	OD	Conc	Factor Range		Slope Check			
				Low	High				
Point 1:	#		0.0	-999999	999999	Allowable Range Check			
Point 2:	#		100.0 £	-999999	999999	<input type="checkbox"/> Reagent Blank			
Point 3:						<input type="checkbox"/> Calibration			
Point 4:						Advanced Calibration			
Point 5:						Operation			
Point 6:						Interval (RB/ACAL)			
Point 7:						No			
Point 8:									
Point 9:									
Point10:									
<Point Cal. For Master Curve>		No. of Correction Points		Use Master Curve		<input type="checkbox"/> Lot Calibration			
	Calibrator	OD	Conc	OD Range		Stability			
				Low	High	Reagent Blanks			
Point 1:						Calibration			
Point 2:						Day			
MB Type Factor:			1-Point Calibration Point			<input type="checkbox"/> With CONC-0			

User Defined

£ For conversion to g/L change the calibrator concentration to 1 and the units to g/L

DRI Ethyl Alcohol Assay - Quantitative

Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Tests	Range
Test Name: # ▾		< >		Type: # ▾	Operation: Yes ▾
Sample Volume	Volume	9.0	μL	Dilution	0 ▾ μL
	Pre-Dilution Rate	1	▾		OD Limit
					Min OD -2.00 Max OD 3.00
Reagent Volume	R1(R1-1)	75	μL	Dilution	0 μL
	R1-2		μL	Dilution	
	R2(R2-1)	75	μL	Dilution	0 μL
					Reagent OD Limit
					First Low -2.00 High 3.00
					Last Low -2.00 High 3.00
Common Reagent	Type	None		Name	
Wavelength:	Pri.	340	▾nm	Sec.	410 ▾nm
Method:		FIXED1*	▾		
Reaction slope:		+	▾	Dynamic Range	Low 10 High 600
Measuring Point 1	First	13		Last	17
				Correlation Factor	A 1 B 0
Measuring Point 2	First			Last	
				Factor for Maker	A 1 B 0
Linearity:			%		
Lag Time Check		No	▾	On-board Stability Period	# Day # Hour

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Tests	Range
Test Name: # ▾		< >		Type: # ▾	
Value/Flag: # ▾					
Level		Low	High		
		#	#		
Specific Ranges:					
		From	To	Low	High
	Sex	Year	Month	Year	Month
□ 1.	# ▾	#	#	#	#
□ 2.	# ▾	#	#	#	#
□ 3.	# ▾	#	#	#	#
□ 4.	# ▾	#	#	#	#
□ 5.	# ▾	#	#	#	#
□ 6.	# ▾	#	#	#	#
7. Standard Demographics					
8. Not within expected values					
Panic Value	Low	#	High	#	Unit mg/dL
					Decimal Places #

*Can be run as RATE1
User Defined Values

DRI Ethyl Alcohol Assay - Quantitative
Beckman Coulter System Parameters, AU5800, continued

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

User Defined Values

£ For conversion to g/L change the calibrator concentration to 1 and the units to g/L

© 2017 Thermo Fisher Scientific, Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AU Series Systems are the registered trademarks of Beckman Coulter.

End