

DRI[®] CREATININE-DETECT ASSAY APPLICATION
BECKMAN COULTER AU5800[®]

Catalog No. 1797

Intended for the quantitative determination of creatinine in human urine for the detection of adulteration by dilution or substitution with a non-urine solution.

For In Vitro Diagnostic and Rx Use

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

Item	Size	Catalog Number
DRI [®] Creatinine-Detect Assay	500 mL	1797
DRI [®] Creatinine-Detect Calibrator Kit	2 x 25 mL	100272
DRI [®] Creatinine-Detect 1.3 mg/dL Control	25 mL	100273
DRI [®] Creatinine-Detect 7.5 mg/dL Control	25 mL	100274
DRI [®] Creatinine-Detect 23.0 mg/dL Control	25 mL	100275

To place an order or for technical service, contact:

USA	In Europe
Tel: (800) 232-3342 Fax: (510) 829-8115	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) 829-8115 / Fax: (510) 979-5420

B·R·A·H·M·S GmbH, Neuendorfstrasse 25, 16761, Hennigsdorf, Germany
Tel: +49 (0)851-88 6890/ Fax: +49 (0)851-88 68910

Continued on next page

**Reagent
Storage**

Refer to the package insert for information on reagent storage.

**Procedure for
Analyzer**

1. Refer to the operator's manuals for information on analyzer operation.
 2. See package insert for specimen preparation.
 3. Dispense adequate amounts of Reagent 1 and Reagent 2 into appropriate containers.
 4. Place the filled reagent containers on the reagent trays in positions defined by the user. **Ensure that reagents have equilibrated to temperature of analyzer reagent compartment before starting analysis.**
 5. Perform a reagent volume check as instructed in the operator's manual.
 6. Select "All" to calibrate the test. It is not necessary to test the blank rack during calibration.
 7. Define a worklist as instructed in the operator's manual. Press START to begin analysis.
 8. Results for calibrators will be printed in Optical Density (OD). Results for samples will be printed in mg/dL.
-

**Results and
Data
Interpretation**

Results for samples will be printed in mg/dL.

**Beckman Coulter AU5800 Parameters
DRI Creatinine-Detect Assay**

Parameters		Specific Test Parameters					
General	LIH	ISE	HbA1c		Calculated Test	Range	
Test Name:		#	<	>	Type: Urine	Operation: Yes	
Sample Volume	2.5	μL	Dilution	0	μL	OD Limit	
Pre-Dilution Rate	1	▽	Diluent Bottle	#	▽	Min.OD: -2.00 Max.OD: 3.00	
Rgt. Volume	R1(R1-1)	80	μL	Dilution	0	μL	
	R1-2		μL	Dilution		μL	
	R2(R2-1)	80	μL	Dilution	0	μL	
Common Rgt. Type	None		Name	None		Dynamic Range Low: 0.1 High: 184	
Wavelength	Pri	520	▽nm	Sec.	600	▽nm	
Method	FIXED1*		Correlation Factor A	1		High B: 0	
Reaction Slope	+		Factor for Maker A	1		High B: 0	
Measuring Point1 1 st	13		Last	17		Onboard Stability Period: # Day # Hour	
Measuring Point2 1 st			Last			LIH Influence Check: # ▽	
Linearity Limit						Lipemia: ▽	
Lag Time Check	No					Icterus: ▽	
						Hemolysis: ▽	

Parameters		Specific Test Parameters					
General	LIH	ISE	HbA1c		Calculated Tests	Range	
Test Name		#	<	>	Type: Urine	▽	
Value/Flag	#				Low: -9999999	High: 9999999	
Level					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	Standard Demographics		#	#		
	8	Not within expected values		#	#		
Panic Value	Low	#	High	#	Unit: mg/dL	Decimal Places: #	

User defined

Beckman Coulter AU5800 Parameters
DRI Creatinine-Detect Assay (continued)

Parameters		Calibration Parameters			
Calibrators		Calibration Specific		STAT Table Calibration	
Test Name	# ▾	<	>	Type	Urine ▾ <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	# ▾		2.0	-99999	99999
Point-2	# ▾		20.0	-99999	99999
Point-3	▾				
Point-4	▾				
Point-5	▾				
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		▾
	Calibrator	OD	Conc	Low	High
Point-1	▾				
Point-2	▾				
MB Type Factor		1-Point Calibration Point		None ▾ <input type="checkbox"/> with Conc-0	
				Stability	
				Reagent Blank	# Day # Hour
				Calibration	# Day # Hour

User defined

© 2017 Thermo Fisher Scientific Inc. All rights reserved.
All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.
CEDIA is a trademark of Roche Diagnostics.
AU Series Systems are trademarks of Beckman Coulter.

End