

CEDIA® OPIATE APPLICATION
BECKMAN COULTER AU480®, AU680®



Catalog No. 100089, 100098, 1661248, 10016429

Intended for the Qualitative and Semiquantitative Determination of Opiate in Human Urine

For In Vitro Diagnostic Use 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.

Ordering Information Materials available from Microgenics, a part of Thermo Fisher Scientific:

Item	Size	Catalog Number
CEDIA Opiate Assay Reagents	3x17 mL	10016429
	3x17 mL	100089
	65 mL	100098
	495 mL	1661248
CEDIA Negative Calibrator	5 mL	1557416
	15 mL	1661388
CEDIA Primary Calibrator	5 mL	1815326
	15 mL	1815334
CEDIA Secondary Calibrator	5 mL	1730428
	15 mL	1730517
CEDIA Intermediate Calibrator	5 mL	1730380
	15 mL	1732218
CEDIA High Calibrator	5 mL	1730398
	15 mL	1732226
MGC Clinical DAU Control Set	3 x 5 mL	100201

To place an order or 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

Thermo Fisher Scientific Oy, Ratastie 2, P.O. Box 100, 01621 Vantaa, Finland

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 Reagent 1 (EA reagent) and Reagent 2 (ED Reagent) into appropriate containers. **Ensure that reagents have equilibrated to temperature of analyzer reagent compartment before starting analysis.**

Results and Data Interpretation

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

CEDIA Opiate Assay – Qualitative (300 ng/mL cutoff) Beckman Coulter AU480 & AU680 Parameters

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	#	<	>	Type:	Urine	Operation:	Yes			
Sample Volume	2.0	μL	Dilution	0	μL	OD Limit				
Pre-Dilution Rate	1		Min. OD	-2.00	Max.	3.00				
Reagents	R1(R1-	87	μL	Dilution	0	μL	Reagent OD			
							First Low	-2.00	High	3.00
							Last Low	-2.00	High	3.00
	R2 (R2-1)	87	μL	Dilution	0	μL	Dynamic Range	#	High	#
							Correlation Factor A	1	B	†
							Factor for Maker A	1	B	0
Wavelength:	Pri.	570	nm	Sec.	660	nm	Onboard Stability	#	Days	#
Method:		FIXED*					LIH Influence Check	#		
Reaction slope:		+					Lipemia			
Measuring Point 1:	First	24		Last	27		Icterus			
Measuring Point 2:	First			Last			Hemolysis			
Linearity:			%							
No Lag Time:		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	Low	High		
	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	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 2.0 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.

**CEDIA Opiate Assay – Qualitative (300 ng/mL cutoff)
Beckman Coulter AU480 & AU680 Parameters, *continued***

(Option 1)

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

(Option 2 or 3)

Parameters	Calibration Parameters		
Calibrators	Calibration Specific	STAT Table Calibration	
Test Name <input type="text" value="#"/> ▾	<input type="text" value="<"/> <input type="text" value=">"/>	Type <input type="text" value="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 >		Factor Range	Slope Check <input type="text" value="+"/> ▾
	Calibrator	OD	Conc
			Low High
Point-1	<input type="text" value="#"/>	<input type="text"/>	<input type="text" value="100"/>
Point-2	<input type="text"/>	<input type="text"/>	<input type="text" value="-9999999"/>
Point-3	<input type="text"/>	<input type="text"/>	<input type="text" value="9999999"/>
Point-4	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-5	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-6	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-7	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-8	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-9	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-10	<input type="text"/>	<input type="text"/>	<input type="text"/>
< Point Cal. For Master Curve >		No. of Correction Points <input type="text"/>	Use Master Curve <input type="text"/>
	Calibrator	OD	Conc
			Low High
Point-1	<input type="text"/>	<input type="text"/>	<input type="text"/>
Point-2	<input type="text"/>	<input type="text"/>	<input type="text"/>
MB Type Factor <input type="text"/>		1-Point Calibration Point <input type="text"/>	<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 <input type="text" value="No"/> ▾
		Interval (RB/ACAL)	<input type="text"/>
		<input type="checkbox"/> Lot Calibration	
		Allowable Range Check	<input type="checkbox"/> Reagent Blank <input type="text"/>
			<input type="checkbox"/> Calibration <input type="text"/>

CEDIA Opiate Assay – Semiquantitative (300 ng/mL cutoff) Beckman Coulter AU480 & AU680 Parameters

Specific Test Parameters										
General		LIH		ISE		Range				
Test Name:	#	<	>	Type:	Urine	Operation:	Yes			
Sample Volume	2.0	μL	Dilution	0	μL	OD Limit				
Pre-Dilution Rate	1					Min. OD	-2.00	Max.	3.00	
Reagents	R1(R1-	87	μL	Dilution	0	μL	Reagent OD			
						First Low	-2.00	High	3.00	
						Last Low	-2.00	High	3.00	
	R2 (R2-1)	87	μL	Dilution	0	μL	Dynamic Range	#	High	#
Wavelength:	Pri.	570	nm	Sec.	660	nm	Correlation Factor A	1	B	0
Method:	FIXED1*									
Reaction slope:	+									
Measuring Point 1:	First	24	Last	27	LIH Influence Check	#				
Measuring Point 2:	First		Last		Lipemia					
Linearity:										
No Lag Time:	No									
					Icterus					
					Hemolysis					

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

