

CEDIA[®] AMPHETAMINE OFT ASSAY APPLICATION (ANZ) BECKMAN COULTER AU680[®]

Catalog No. 10011931, 10011932

The Thermo Scientific CEDIA Amphetamine OFT Assay is intended for use in the semiquantitative determination of amphetamine in human oral fluid at a cutoff concentration of 17 ng/mL (diluted oral fluid) or 51 ng/mL (neat oral fluid). The specimen must be collected exclusively with the Oral-Eze[®] Oral Fluid Collection System.

For In Vitro Diagnostic Use Only
For International Use Only – Not for sale in US
(For Use with Oral-Eze Oral Fluid Collection System)

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 Amphetamine OFT Assay	3 x 18 mL 1 x 65 mL	10011931 10011932
CEDIA Multi-Drug OFT Negative Calibrator	1 x 20 mL	10016864
CEDIA Multi-Drug OFT Calibrator 1	1 x 10 mL	10016882
CEDIA Multi-Drug OFT Calibrator 2	1 x 10 mL	10016883
CEDIA Multi-Drug OFT Calibrator 3	1 x 10 mL	10016884
CEDIA Multi-Drug OFT Control 1	1 x 15 mL	10017711
CEDIA Multi-Drug OFT Control 2	1 x 15 mL	10017712
CEDIA Multi-Drug OFT Control 3	1 x 15 mL	10017713

To place an order or for technical service, contact:

In Europe
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 package insert for information on reagent storage.

Analyzer Procedure

1. Set up the Beckman Coulter AU680 as instructed in the operator's manual.
2. Dispense adequate amounts of calibrators/control and sample(s) into sample cups in the appropriate racks.
3. Dispense adequate amounts of EA and ED reagents into appropriate containers.
4. Place the filled reagent containers on the reagent trays in positions defined by the user. **EA Reagent is used as R1; ED Reagent is used as R2. Make sure the reagents have equilibrated to the temperature of the analyzer reagent compartment before starting analysis.**
5. Perform a reagent volume check as instructed in the operator's manual.
6. Define a worklist as instructed in the operator's manual. Press START to begin 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

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

CEDIA Amphetamine OFT Assay (ANZ), Semiquantitative Beckman Coulter System Parameters, AU680

Specific Test Parameters																			
General		LIH		ISE		Range													
Test Name:		AmpOFT ▾		< >		Type:		Urine ▾		Operation:		Yes ▾							
Sample Volume		25 μL		Dilution		0 μL		OD Limit											
Pre-Dilution Rate		1		Min. OD		-2.00		Max. OD				3.00							
Reagents		R1(R1-		75 μL		Dilution		0 μL		Reagent OD									
								First Low		-2.00		High		3.00					
								Last Low		-2.00		High		3.00					
		R2 (R2-1)		75 μL		Dilution		0 μL		Dynamic Range				-9999.9		High		9999.9	
										Correlation Factor A		*		B		0			
Wavelength:		Pri.		570 nm		Sec.		660 nm		Factor for Maker A		1		B		0			
Method:		FIXED1 ▾								Onboard Stability		#		Days		#		Hour	
Reaction slope:		+								LIH Influence Check		#							
Measuring Point 1:		First		24		Last		27		Lipemia				▽					
Measuring Point 2:		First				Last				Icterus				▽					
Linearity:				%						Hemolysis				▽					
No Lag Time:		No																	

For Specific Test Parameters → Range Tab, enter “ng/mL” for Unit and “1” for Decimal Places.

* Option 1: When the correlation factor “1” is used, the sample results represent diluted oral-fluid concentrations.

* Option 2: When the correlation factor “3” is used, the sample results are multiplied by 3 to represent undiluted (neat) oral-fluid concentrations.

Parameters		Specific Test Parameters													
General		LIH		ISE		Calculated Tests				Range					
Test Name		# ▾		< >		Type		Urine ▾							
Value/Flag		# ▾													
Level				Low		High		Panic Value							
				-9999999		‡									
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		ng/mL		Decimal Places		#									

**CEDIA Amphetamine OFT Assay (ANZ), Semiquantitative
Beckman Coulter System Parameters, AU680, *continued***

Parameters		Calibration Parameters					
Calibrators		Calibration Specific		STAT Table Calibration			
Test Name	<input type="text" value="AmpOFT"/> ▾	<input type="text" value="<"/>	<input type="text" value=">"/>	Type	<input type="text" value="Serum"/> ▾ <input type="checkbox"/> Use Serum Cal.		
Calibration Type	<input type="text" value="4AB"/> ▾	Formula	<input type="text" value="Polygonal"/> ▾	Counts	<input type="text" value="2"/> ▾		
< Calibrator Parameters >			Range		Slope Check	<input type="text" value="+"/> ▾	
	Calibrator	OD	Conc	Low	High	Allowable Range Check	
Point-1	<input type="text" value="1."/> ▾		0.0	-2.0000	3.0000	<input type="checkbox"/> Reagent Blank	
Point-2	<input type="text" value="2."/> ▾		10.0	-2.0000	3.0000	<input type="checkbox"/> Calibration	
Point-3	<input type="text" value="3."/> ▾		17.0	-2.0000	3.0000		
Point-4	<input type="text" value="4."/> ▾		40.0	-2.0000	3.0000		
Point-5	<input type="text" value=""/> ▾					Advanced Calibration	
Point-6	<input type="text" value=""/> ▾					Operation	
Point-7	<input type="text" value=""/> ▾					Interval (RB/ACAL)	<input type="text" value="No"/> ▾
Point-8	<input type="text" value=""/> ▾						
Point-9	<input type="text" value=""/> ▾					<input type="checkbox"/> Lot Calibration	
Point-10	<input type="text" value=""/> ▾						
< Point Cal. For Master Curve >		No. of Correction Points	<input type="text" value=""/> ▾	Use Master Curve	<input type="text" value=""/> ▾		
	Calibrator	OD	Conc	OD Range		Stability	
Point-1	<input type="text" value=""/> ▾			Low	High	Reagent Blank	<input type="text" value=""/> Day <input type="text" value=""/> Hour
Point-2	<input type="text" value=""/> ▾					Calibration	<input type="text" value=""/> Day <input type="text" value=""/> Hour
MB Type Factor	<input type="text" value=""/> <input type="text" value=""/>	1-Point Calibration Point	<input type="text" value="None"/> ▾	<input type="checkbox"/> with Conc-0			

© 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