

CEDIA® PHENCYCLIDINE (PCP) APPLICATION CE
**Beckman Coulter AU400®, AU480®, AU640®, AU680®,
AU2700®, AU5400®, AU5800®**

Catalog No. 100172, 100173, 1815784

Intended for the Qualitative and Semiquantitative Determination of Phencyclidine 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.

Materials available from Microgenics, a part of Thermo Fisher Scientific

Ordering Information

Item	Size	Thermo Fisher Scientific Catalog No.
CEDIA Phencyclidine Assay Reagents	3x17 mL	100172
	65 mL	100173
	495 mL	1815784
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 Primary DAU Control Set	3 x 5 mL	100200

To place an order or for technical service contact:

In USA	In Europe
(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

EC REP Thermo Fisher Scientific Oy, Ratastie 2, P.O.Box 100, 01621 Vantaa, Finland
Tel: +358-9-329100/Fax: +358-9-32910300

Continued on next page

Reagent Storage

Refer to the package insert for information on reagent storage.

Procedure for Analyzer

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

Results for samples will be printed in ng/mL

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

Additional, user-definable parameters present results that are interpretable in different ways. For further information, contact Microgenics Technical Support or your local Microgenics representative.

CEDIA Phencyclidine Assay - Qualitative

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

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	PCPQ	<	>	Type:	Urine	Operation:	Yes			
Sample:	Volume	4.0	μL	Dilution	0	μL	Pre-Dilution Rate:	1		
Reagents:	R1 Volume	87	μL	Dilution	0	μL	Min OD	Max OD		
	R2 Volume	87	μL	Dilution	0	μL	L	-2.00	H	2.50
Wavelength:	Pri.	570	▼	Sec.	660	▼	Reagent OD limit:			
Method:	RATE* ▼			First L	-2.00	First H	2.50			
				Last L	-2.00	Last H	2.50			
Reaction slope:	+ ▼			Dynamic Range:						
Measuring Point 1:	First	24	Last	27	L	#	H	#		
Measuring Point 2:	First		Last		Correlation Factor:					
Linearity:				A	1	B	0			
No Lag Time:	No ▼			On-board stability period:	#					

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	PCPQ	<	>	Type:	Urine					
Value/Flag:	#	Level L:	#	Level H:	#					
Normal Ranges:		Age L		Age H		L		H		
<input type="checkbox"/>	1.	Sex	Year	Month	Year	Month	#	#	#	#
<input type="checkbox"/>	2.	#	#	#	#	#	#	#	#	#
<input type="checkbox"/>	3.	#	#	#	#	#	#	#	#	#
<input type="checkbox"/>	4.	#	#	#	#	#	#	#	#	#
<input type="checkbox"/>	5.	#	#	#	#	#	#	#	#	#
<input type="checkbox"/>	6.	#	#	#	#	#	#	#	#	#
7. None Selected						#		#		
8. Out of Range		L		H		#		#		
Panic Value:	#		#		Unit:	#		Decimal places: #		

Calibration Specific										
General		ISE								
Test Name:	PCPQ	<	>	Type:	Urine					
Calibration Type:	AB	Formula:	Y=AX+B	Counts:	2	Process:	CONC			
Point 1:	Cal. No.	OD	CONC	Factor/OD-L	Factor/OD-H					
	#		100	-99999	99999					
Point 2:										
Point 3:										
Point 4:										
Point 5:										
Point 6:										
Point 7:										
1-Point Cal. Point:	<input type="checkbox"/>	<input type="checkbox"/>	With CONC-0	Slope Check	+	Advanced Calibration				
MB Type Factor:			Calibration Stability Period:	#						

User Defined
 *Can also be run as FIXED

CEDIA Phencyclidine Assay - Semiquantitative

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

Specific Test Parameters											
General		LIH	ISE	Range							
Test Name:	PCPSQ ▾			<	>	Type:	Urine ▾		Operation:	Yes ▾	
Sample:	Volume	4.0	μL	Dilution	0	μL	Pre-Dilution Rate:	1			
Reagents:	R1 Volume	87	μL	Dilution	0	μL	Min OD	Max OD			
	R2 Volume	87	μL	Dilution	0	μL	L	-2.00	H 2.50		
Wavelength:	Pri.	570	▾	Sec.	660	▾	Reagent OD limit:	First L	-2.00	First H	2.50
Method:	RATE1* ▾					Last L	-2.00	Last H 2.50			
Reaction slope:	+ ▾					Dynamic Range:	L	0	H 150		
Measuring Point 1:	First	24		Last	27		Correlation Factor:	A	1	B	0
Measuring Point 2:	First			Last			On-board stability period:	#			
Linearity:						%					
No Lag Time:	No					▾					

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	PCPSQ ▾			<	>	Type:	Urine ▾			
Value/Flag:	# ▾		Level L:	#		Level H:	#			
Normal Ranges:	Age L		Age H		L		H			
	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. None Selected				#		#			
	8. Out of Range		L		H		#		#	
Panic Value:	#		#		Unit:	ng/mL		Decimal places: #		

Calibration Specific											
General		ISE									
Test Name:	PCPSQ ▾			<	>	Type:	Urine ▾				
Calibration Type:	4AB ▾		Formula:	POLYGONAL ▾		Counts:	2		Process:	CONC ▾	
Point 1:	Cal. No.	#	OD		CONC	0	Factor/OD-L	-2.00	Factor/OD-H		2.50
Point 2:	Cal. No.	#			25		-2.00	2.50			
Point 3:	Cal. No.	#			75		-2.00	2.50			
Point 4:	Cal. No.	#			150		-2.00	2.50			
Point 5:	Cal. No.										
Point 6:	Cal. No.										
Point 7:	Cal. No.										
1-Point Cal. Point:	<input type="checkbox"/>	<input type="checkbox"/>	With CONC-0	Slope Check	+	▾	Advanced Calibration	# ▾			
MB Type Factor:						Calibration Stability Period:	#				

*Can also be run as FIXED1
#User Defined

CEDIA Phencyclidine Assay - Qualitative

Beckman Coulter System Parameters, AU480/AU680

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	PCPQ ▾	<	>	Type	Urine ▾	Operation	Yes ▾		
Sample Volume	4.0 μL	Dilution	0 ▾ μL	OD Limit		Min.OD	-2.0000	Max.OD	3.0000
Pre-Dilution Rate	1 ▾								
Reagent Volume	R1(R1-1) 87 μL	Dilution	0 μL	Reagent OD Limit		First	Low -2.0000	High	3.0000
	R2(R2-1) 87 μL	Dilution	0 μL			Last	Low -2.0000	High	3.0000
Common Reagent	Type None	Name		Dynamic Range		Low	#	High	#
Wavelength	Pri. 570 ▾ nm	Sec.	660 ▾ nm	Correlation Factor		A	1	B	0
Method	FIXED* ▾			Factor for Maker		A	1	B	0
Reaction slope	+ ▾								
Measuring Point-1	First 24	Last	27	Onboard Stability Period			30 Day	0 Hour	
Measuring Point-2	First	Last							
Linearity Limit									
Lag Time Check	No ▾								

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c	Calculated Tests	Range				
Test Name	PCPQ ▾	<	>	Type	Urine ▾				
Value/Flag	# ▾								
Level		Low	#	High	#	Panic Value			
						Low	#	High	#
Specific Ranges:		From		To					
	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	#						

CEDIA Phencyclidine Assay - Qualitative

Beckman Coulter System Parameters, AU480/AU680, *continued*

Parameters		Calibration Parameters			
Calibrators		Calibration Specific		STAT Table Calibration	
Test Name	<input type="text" value="PCPQ"/> ▾	<input type="button" value=" <"/>	<input type="button" 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
	Calibrator	OD	Conc	Low	High
Point-1	# ▾		100	-9999999	9999999
Point-2	▾				
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	<input type="text" value="1"/> ▾	Use Master Curve <input type="text" value="1"/> ▾	
			OD Range		
	Calibrator	OD	Conc	Low	High
Point-1	▾				
Point-2	▾				
MB Type Factor		<input type="text"/>	1-Point Calibration Point	<input type="text" value="1"/> ▾	<input type="checkbox"/> with Conc-0
				Stability	
				Reagent Blank	<input type="text" value="30"/> Day <input type="text" value="0"/> Hour
				Calibration	<input type="text" value="0"/> Day <input type="text" value="0"/> Hour
				Advanced Calibration	<input type="text" value="No"/> ▾
				Operation	<input type="text" value="1"/> ▾
				Interval (RB/ACAL)	<input type="text" value="1"/> ▾
				<input type="checkbox"/> Lot Calibration	
				Allowable Range Check	<input type="text" value="None"/> ▾
				<input type="checkbox"/> Reagent Blank	<input type="text"/>
				<input type="checkbox"/> Calibration	<input type="text"/>

User defined
 * Can also be run as RATE

CEDIA Phencyclidine Assay - Semiquantitative Beckman Coulter System Parameters, AU480/AU680

Specific Test Parameters										
General		LIH	ISE	Range						
Test Name:	PCPSQ	<	>	Type:	Urine	Operation:	Yes			
Sample Volume	4.0	μL	Dilution	0	μL	OD Limit				
Pre-Dilution Rate	1					Min. OD	-2.00	Max. OD	3.00	
Reagents Volume:	R1(R1-1)	87	μL	Dilution	0	μL	Reagent OD limit:			
						First Low	-2.00	High	3.00	
						Last Low	-2.00	High	3.00	
	R2 (R2-1)	87	μL	Dilution	0	μL	Dynamic Range Low	0	High	150
						Correlation Factor A	1	B	0	
Wavelength:	Pri.	570	nm	Sec.	660	nm	Factor for Maker 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	▾			

Specific Test Parameters									
General		ISE	Range						
Test Name:	PCPSQ	<	>	Type:	Urine				
Value/Flag:	VALUE	Level L:	#	Level H:	#				
Specific Ranges:									
	Sex	Year	Month	Year	Month	Low	High	Panic Value	
<input type="checkbox"/>	1.	#	#	#	#	#	#	Low	High
<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	#					

User defined
* Can also be run as RATE1

CEDIA Phencyclidine Assay - Semiquantitative Beckman Coulter System Parameters, AU480/AU680, *continued*

Calibration Specific										
General		ISE								
Test Name:		PCPSQ ▾			< >		Type		Urine ▾	<input type="checkbox"/> Use Serum Cal.
Calibration Type:		4AB ▾	Formula:		Polygonal ▾		Counts:		2	
<Calibrator Parameters>										
	Calibrator †	OD	Conc	Factor Range			Slope Check		+ ▾	
				Low	High		Allowable Range Check			
Point 1:	# ▾		0	-2.00	3.00		<input type="checkbox"/> Reagent Blank			
Point 2:	# ▾		25	-2.00	3.00		<input type="checkbox"/> Calibration			
Point 3:	# ▾		75	-2.00	3.00		Advanced Calibration			
Point 4:	# ▾		150	-2.00	3.00		Operation		No ▾	
Point 5:							Interval (RB/ACAL)			
Point 6:										
Point 7:										
Point 8:										
Point 9:										
Point10:										
<Point Cal. For Master Curve>										
	Calibrator	No. of Correction Points		Use Master Curve			<input type="checkbox"/> Lot Calibration			
		OD	Conc	OD Range			Stability			
				Low	High		Reagent Blanks Calibration		# Day # Hour	
Point 1:							#		Day	#
Point 2:							30		Day	0
	MB Type Factor:		1-Point Calibration Point	None ▾			<input type="checkbox"/> With CONC-0			

User defined

CEDIA Phencyclidine Assay – Qualitative

Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters					
General	LIH	ISE	HbA1c		Calculated Test	Range	
PCPSQ ▾		< >		Type: Urine ▾	Operation Yes ▾		
Sample Volume	4 μL	Dilution	0 μL	OD Limit			
Pre-Dilution Rate	1 ▾	Diluent Bottle	Outside ▾	Min.OD	-2.00	Max.OD	3.00
Rgt. Volume	R1(R1-1) 87 μL	Dilution	0 μL	Reagent OD Limit			
	R1-2	Dilution		1 st .	Low -2.00	High 3.00	
	R2(R2-1) 87 μL	Dilution	0 μL	Last	Low -2.00	High 3.00	
Common Rgt. Type	None	Name	None	Dynamic Range Low	#	High	#
Wavelength	Pri 570 ▾nm	Sec.	660 ▾nm	Correlation Factor A	1	B	0
Method	FIXED* ▾			Factor for Maker A	1	B	0
Reaction Slope	+ ▾			Onboard Stability Period	30 Day	0 Hour	
Measuring Point1 1 st	24	Last	27	LIH Influence Check	# ▾		
Measuring Point2 1 st		Last		Lipemia		▾	
Linearity Limit	%			Icterus		▾	
Lag Time Check	No ▾			Hemolysis		▾	

Parameters		Specific Test Parameters					
General	LIH	ISE	HbA1c		Calculated Test	Range	
Test Name: PCPQ ▾		< >		Type: Urine ▾			
Value/Flag: # ▾				Level	Low #	High #	
Specific Ranges:		From		To	Low	High	
<input type="checkbox"/> 1.	Sex # ▾	Year #	Month #	Year #	Month #	Low #	High #
<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 #	Decimal Places #			

User defined.
 * Can also be run as RATE

CEDIA Phencyclidine Assay – Qualitative Beckman Coulter System Parameters, AU5800, *continued*

Parameters		Calibration Parameters	
Calibrators	Calibration Specific	STAT Table Calibration	
Test Name: <input type="text" value="PCPQ"/> ▾		Type <input type="text" value="Urine"/> ▾	
<input type="checkbox"/> Use Serum Cal		Cuvette <input type="text" value=""/> ▾	
Calibration Type: <input type="text" value="AB"/> ▾		Formula: <input type="text" value="Y=AX+B"/> ▾	
< Calibrator Parameters >		Counts: <input type="text" value="2"/> ▾	
Calibrator	OD	Conc	OD Range
			Low High
Point-1	# ▾	100	-99999 99999
Point-2	▾		
Point-3	▾		
Point-4	▾		
Point-5	▾		
Point-6	▾		
Point-7	▾		
Point-8	▾		
Point-9	▾		
Point-10	▾		
		Slope Check <input type="text" value="+"/> ▾	
		Allowable Range Check	
		<input type="checkbox"/> Reagent Blank <input type="text" value=""/>	
		<input type="checkbox"/> Calibration <input type="text" value=""/>	
		Advanced Calibration	
		Operation <input type="text" value="No"/> ▾	
		Interval (RB/ACAL) <input type="text" value=""/> ▾	
		<input type="checkbox"/> Lot Calibration	
<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
			Low High
Point-1	▾		
Point-2	▾		
MB Type Factor <input type="text" value=""/>		1-Point Calibration Point <input type="text" value="None"/> ▾	
		<input type="checkbox"/> with Conc-0	
		Stability	
		Reagent Blank <input type="text" value="#"/> Day <input type="text" value="#"/> Hour	
		Calibration <input type="text" value="30"/> Day <input type="text" value="0"/> Hour	

User Defined Values

CEDIA Phencyclidine Assay – Semiquantitative Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: <input type="text" value="PCPSQ"/> <input type="button" value="<"/> <input type="button" value=">"/>		Type: <input type="text" value="Urine"/>		Operation <input type="text" value="Yes"/>	
Sample Volume	<input type="text" value="4"/> μL	Dilution	<input type="text" value="0"/> μL	OD Limit	
Pre-Dilution Rate	<input type="text" value="1"/>	Diluent Bottle	<input type="text" value="#"/>	Min.OD	<input type="text" value="-2.00"/> Max.OD <input type="text" value="3.00"/>
Rgt. Volume	R1(R1-1) <input type="text" value="87"/> μL	Dilution	<input type="text" value="0"/> μL	Reagent OD Limit	
	R1-2 <input type="text" value=""/>	Dilution	<input type="text" value=""/>	1 st . Low	<input type="text" value="-2.00"/> High <input type="text" value="3.00"/>
	R2(R2-1) <input type="text" value="87"/> μL	Dilution	<input type="text" value="0"/> μL	Last Low	<input type="text" value="-2.00"/> High <input type="text" value="3.00"/>
Common Rgt. Type	<input type="text" value="None"/>	Name	<input type="text" value="None"/>	Dynamic Range Low	<input type="text" value="0"/> High <input type="text" value="150"/>
Wavelength	Pri <input type="text" value="570"/> nm	Sec.	<input type="text" value="660"/> nm	Correlation Factor A	<input type="text" value="1"/> B <input type="text" value="0"/>
Method	<input type="text" value="FIXED1*"/>	Factor for Maker A		<input type="text" value="1"/> B <input type="text" value="0"/>	
Reaction Slope	<input type="text" value="+"/> <input type="button" value="v"/>	Onboard Stability Period		<input type="text" value="30"/> Day	<input type="text" value="0"/> Hour
Measuring Point1 1 st	<input type="text" value="24"/>	Last	<input type="text" value="27"/>	LIH Influence Check <input type="text" value="#"/>	
Measuring Point2 1 st	<input type="text" value=""/>	Last	<input type="text" value=""/>	Lipemia <input type="text" value="v"/>	
Linearity Limit	<input type="text" value=""/>			Icterus <input type="text" value="v"/>	
Lag Time Check	<input type="text" value="No"/>			Hemolysis <input type="text" value="v"/>	

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: <input type="text" value="PCPSQ"/> <input type="button" value="<"/> <input type="button" value=">"/>		Type: <input type="text" value="Urine"/>			
Value/Flag: <input type="text" value="#"/>		Level To		Low <input type="text" value="#"/>	High <input type="text" value="#"/>
Specific Ranges:		From	To	Low	High
	Sex	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="text" value="#"/>
<input type="checkbox"/> 2.	<input type="text" value="#"/>	<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="text" value="#"/>
<input type="checkbox"/> 4.	<input type="text" value="#"/>	<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="text" value="#"/>
<input type="checkbox"/> 6.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>
	7. Standard demographics			<input type="text" value="#"/>	<input type="text" value="#"/>
	8. Not within expected values			<input type="text" value="#"/>	<input type="text" value="#"/>
Panic Value	Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>	Unit <input type="text" value="ng/mL"/> Decimal Places <input type="text" value="#"/>

* Can be run as RATE1
User Defined Values

CEDIA Phencyclidine Assay – Semiquantitative
Beckman Coulter System Parameters, AU5800, *continued*

Parameters		Calibration Parameters	
Calibrators	Calibration Specific	STAT Table Calibration	

Test Name: < > Type Cuvette
 Use Serum Cal

Calibration Type: Formula: Counts:

< Calibrator Parameters >

	Calibrator	OD	Conc	OD Range		Slope Check
				Low	High	
Point-1	<input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="0"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	<input type="text" value="+"/>
Point-2	<input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="25"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Allowable Range Check <input type="checkbox"/> Reagent Blank <input type="text" value=""/> <input type="checkbox"/> Calibration <input type="text" value=""/>
Point-3	<input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="75"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	
Point-4	<input type="text" value="#"/>	<input type="text" value=""/>	<input type="text" value="150"/>	<input type="text" value="-2.0"/>	<input type="text" value="3.0"/>	Advanced Calibration Operation <input type="text" value="No"/> Interval (RB/ACAL) <input type="text" value=""/> <input type="checkbox"/> Lot Calibration
Point-5	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point-6	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point-7	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point-8	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point-9	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	
Point-10	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	

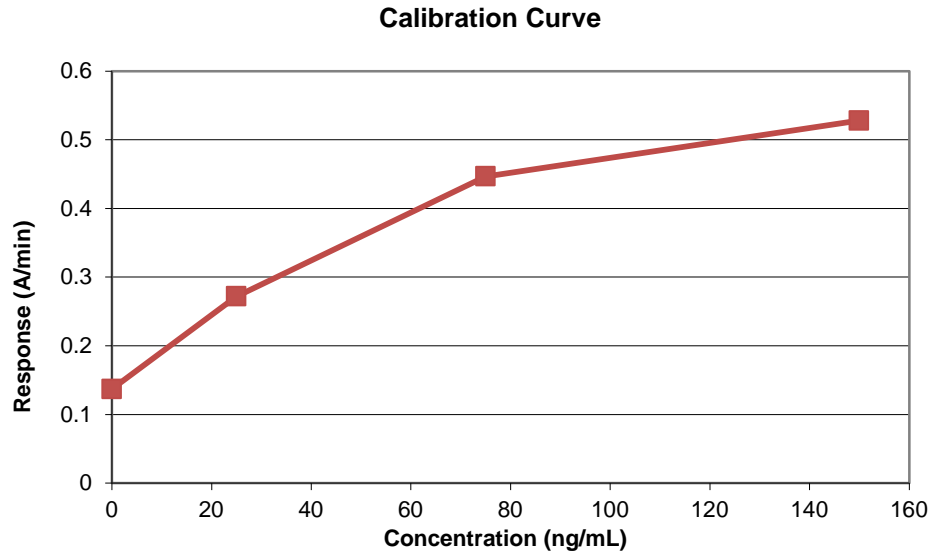
<Point Cal. For Master Curve> No. of Correction Points Use Master Curve

	Calibrator	OD	Conc.	OD Range		Stability
				Low	High	
Point-1	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	Reagent Blank <input type="text" value="#"/> Day <input type="text" value="#"/> Hour Calibration <input type="text" value="30"/> Day <input type="text" value="0"/> Hour
Point-2	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	<input type="text" value=""/>	

MB Type Factor 1-Point Calibration Point with Conc-0

User Defined Values

**Example
Calibration
Curve**



Limit Of Blank The negative calibrator was run against the same calibration curve for 21 replicates. The LOB is calculated as $3 \times SD$.

The observed LDD for Phencyclidine on:

AU480: 0.49 ng/mL
AU680: 0.64 ng/mL
AU5800: 0.92 ng/mL

Precision

Samples of the Low Control, Cutoff Calibrator, High Control were tested in replicates of 2, twice per day for 20 days, total N = 80/Level. The results are presented in the following tables:

	Control Low	Cutoff	Control High
Qualitative AU400*			
Mean (mA/min)	275	321	361
Within-Run SD (mA/min)	3.2	3.4	5.3
Within-Run CV (%)	1.2	1.1	1.5
Total SD (A/min)	7.6	8.3	9.5
Total CV (%)	2.8	2.6	2.6
Qualitative AU480			
Mean (mA/min)	221	263	308
Within-Run SD (mA/min)	1.4	1.8	2.5
Within-Run CV (%)	0.6	0.7	0.8
Total SD (A/min)	7.3	8.7	10.4
Total CV (%)	3.3	3.3	3.4
Qualitative AU640*			
Mean (mA/min)	259	306	345
Within-Run SD (mA/min)	2.8	3.4	3.5
Within-Run CV (%)	1.1	1.1	1.0
Total SD (A/min)	5.0	6.8	7.0
Total CV (%)	1.9	2.2	2.0
Qualitative AU680			
Mean (mA/min)	221	262	307
Within-Run SD (mA/min)	1.7	2.0	2.7
Within-Run CV (%)	0.8	0.7	0.9
Total SD (A/min)	4.2	4.5	5.1
Total CV (%)	1.9	1.7	1.7
Qualitative AU5800			
Mean (mA/min)	220	261	307
Within-Run SD (mA/min)	2.9	4.0	3.9
Within-Run CV (%)	1.3	1.5	1.3
Total SD (A/min)	5.7	7.3	7.8
Total CV (%)	2.6	2.8	2.5

*N=60/Level

**Precision
(Continued)**

	Control Low	Cutoff	Control High
Semiquantitative AU400*			
Mean (ng/mL)	17	23	30
Within-Run SD (ng/mL)	0.4	0.5	0.7
Within-Run CV (%)	2.5	2.0	2.3
Total SD (ng/mL)	0.8	1.1	1.5
Total CV (%)	4.7	4.5	5.1
Semiquantitative AU480			
Mean (ng/mL)	17	25	38
Within-Run SD (ng/mL)	0.3	0.5	0.7
Within-Run CV (%)	1.5	1.9	1.9
Total SD (ng/mL)	0.3	0.6	0.8
Total CV (%)	1.7	2.3	2.0
Semiquantitative AU640*			
Mean (ng/mL)	19	26	33
Within-Run SD (ng/mL)	0.5	0.6	0.8
Within-Run CV (%)	2.4	2.4	2.4
Total SD (ng/mL)	0.7	0.9	1.3
Total CV (%)	3.7	3.5	3.9
Semiquantitative AU680			
Mean (ng/mL)	17	25	38
Within-Run SD (ng/mL)	0.3	0.4	0.8
Within-Run CV (%)	2.0	1.7	2.1
Total SD (ng/mL)	0.5	0.6	1.0
Total CV (%)	2.8	2.6	2.6
Semiquantitative AU5800			
Mean (ng/mL)	17	25	38
Within-Run SD (ng/mL)	0.5	0.8	0.9
Within-Run CV (%)	2.8	3.1	2.3
Total SD (ng/mL)	0.5	1.0	1.1
Total CV (%)	3.1	3.9	2.8

*N=60/Level

**Accuracy and
Correlation**

Patient samples were assayed using the CEDIA Phencyclidine Assay on the Beckman Coulter and tested against the reference analyzer, the Hitachi 717.

Qualitative Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU400	100.0	98.7	99.4
AU480	100.0	98.4	99.2
AU640	100.0	100.0	100.0
AU680	100.0	96.9	98.3
AU5800	100.0	98.4	99.2

**Accuracy and Correlation
(Continued)**

CEDIA Phencyclidine Assay Qualitative Method Comparison

	Hitachi 717 +			Hitachi 717 +			Hitachi 717 +				
		-			-			-			
AU400	+	54	1	AU480	+	57	1	AU640	+	54	0
	-	0	77		-	0	62		-	0	78
	Hitachi 717 +			Hitachi 717 +			Hitachi 717 +				
		-			-			-			
AU680	+	55	2	AU5800	+	56	1				
	-	0	63		-	0	63				

Semiquantitative Method Comparison

	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU400	100.0	98.7	99.4
AU480	98.2	98.4	98.3
AU640	100.0	100.0	100.0
AU680	100.0	100.0	100.0
AU5800	98.1	100.0	99.2

CEDIA Phencyclidine Assay Semiquantitative Method Comparison

	Hitachi 717 +			Hitachi 717 +			Hitachi 717 +				
		-			-			-			
AU400	+	54	1	AU480	+	56	1	AU640	+	54	0
	-	0	77		-	1	62		-	0	78
	Hitachi 717 +			Hitachi 717 +			Hitachi 717 +				
		-			-			-			
AU680	+	57	0	AU5800	+	53	0				
	-	0	63		-	1	66				

© 2016 Thermo Fisher Scientific, Inc. All rights reserved.
 AU Series Systems are the registered trademarks of Beckman Coulter Inc.
 All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries.

End