



DRI[®] Ecstasy Assay Application – BECKMAN COULTER[®] UniCel[®] DxC and Synchron[®] System(s)

Beckman Coulter Reorder Number A39935

Homogeneous enzyme immunoassay for the qualitative or semi-quantitative determination of ecstasy drugs 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 storage, quality control, and additional performance data.**

Ordering Information

Materials available from your local Beckman Coulter Representative:

Item	Beckman Coulter Reorder Number
DRI [®] Ecstasy Assay Kit (100 mL)	A39935
DRI [®] Negative Calibrator	A44121
DRI [®] Ecstasy 250 ng/mL Calibrator	A39936
DRI [®] Ecstasy 500 ng/mL Calibrator	A39937
DRI [®] Ecstasy 750 ng/mL Calibrator	A39938
DRI [®] Ecstasy 1000 ng/mL Calibrator	A40566
MGC Select DAU Control Set	A40567
User-Defined Reagent Cartridge (pkg. of 12)	442835

For Technical Support please contact your local Beckman Coulter Representative.

DRI[®] is a registered trademark of Thermo Fisher Scientific Corporation.

Reagent Storage

Refer to the package insert for information on reagent storage.

NOTE:

It is not recommended to leave the reagent on-board for more than 60 days.

Procedure for Analyzer

Refer to the operator's manuals for information on analyzer operation. Dispense adequate amounts of Antibody/Substrate Reagent A and Enzyme Conjugate Reagent E into appropriate compartments of a User Defined Cartridge (PN 442835) as shown in the table below:

	User Defined Cartridge	
DRI[®] Ecstasy Assay Kit	Compartment A	Compartment B
Antibody/Substrate Reagent A	48 mL	
Enzyme Conjugate Reagent E		18 mL

Calibrator Information

Refer to the package insert for information on calibration.

To use qualitative mode monitoring, calibration shall be carried out every 14 days or as indicated by control recovery.

To use semi-quantitative mode monitoring, calibration shall be carried out every 14 days or as indicated by control recovery.

Application Parameter

Parameters DRI Ecstasy Assay chemistry parameter for **qualitative** mode on Synchron System(s).

INSTRUMENT PARAMETERS: SYNCHRON CX

**UniCel DxC
Synchron LX**

Chemistry Name	Ecstasy	NA ^a
Test Name	XTCX	XTCX
Reaction Type:	Rate 1	Rate 1
Units	ng/mL	mA/min
Decimal Precision/Precision:	X.X	X.X
Reaction Direction:	Positive	Positive
Calculation Factor:	0	1000
Math Model:	Linear	DAT
Cal. Time Limit:	336	336
Number of Calibrators:	2	3
#1	0.0	0.0
#2	500.0	500.0
#3		1000.0
#4		
#5		
#6		
Primary Wavelength:	340 nm	340 nm
Secondary Wavelength:	650 nm	650 nm
Sample Volume:	15 µL	10 µL
REAGENTS:		
Primary Inject (first)/First Inject		
Compartment/Component:	A	A
Volume/Dispense Volume:	200 µL	125 µL
Add Time/Inject Time:	NA ^a	NA ^a
Primary Inject (first)/Second Inject		
Compartment/Component:	None	None
Volume/Dispense Volume:	0 µL	NA ^a
Add Time/Inject Time:	NA ^a	-180 ^{a&b}
Second Inject/Third Inject		
Compartment/Component:	B	B
Volume/Dispense Volume:	75 µL	75 µL
Add Time/Inject Time:	368 sec	276 sec
REAGENTS:		
Blank		
Start Read:	237 sec	292 sec

Continued on next page

INSTRUMENT PARAMETERS:	SYNCHRON CX	UniCel DxC Synchron LX
End Read:	300 sec	308 sec
Reaction 1		
Start Read:	96 sec	336 sec
End Read:	144 sec	396 sec
Reaction 2		
Start Read:	NA ^a	NA ^a
End Read:	NA ^a	NA ^a
USABLE RESULT RANGE:		
Lower Limit:	0.000 ^b	0.000 ^b
Upper Limit:	99999.999 ^b	99999.999 ^b
ERROR DETECTION LIMITS:		
Reagent Blank/Blank		
ABS Low Limit:	-1.5 ^b	-1.5 ^b
ABS High Limit:	1.5 ^b	2.2 ^b
Rate Low Limit:	NA ^a	-1.5 ^b
Rate High Limit:	NA ^a	2.200 ^b
Mean Deviation:	NA ^a	2.200 ^b
Reaction/Reaction 1		
ABS Low Limit:	-1.5 ^b	-1.5 ^b
ABS High Limit:	1.5 ^b	2.2 ^b
Rate Low Limit:	NA ^a	-1.5 ^b
Rate High Limit:	NA ^a	2.200 ^b
Mean Deviation:	NA ^a	2.200 ^b
Reaction 2		
ABS Low Limit:	NA ^a	-1.5 ^b
ABS High Limit:	NA ^a	2.2 ^b
Rate Low Limit:	NA ^a	-1.5 ^b
Rate High Limit:	NA ^a	2.200 ^b
Mean Deviation:	NA ^a	2.200 ^b
SUBSTRATE DEPLETION		
Initial Rate:	99.999	99.999
Delta ABS:	1.5 ^b	2.2 ^b
MULTI POINT SPAN:	0.000 ^b	0.000 ^b

^a NA=Not applicable

^b Denotes default value

Continued on next page

Parameters DRI Ecstasy Assay chemistry parameter for **semi-quantitative** mode on the Synchron System(s).

INSTRUMENT PARAMETERS: SYNCHRON CX

**UniCel DxC
Synchron LX**

Chemistry Name	Ecstasy	NA ^a
Test Name	XTCX	XTCX
Reaction Type:	Rate 1	Rate 1
Units	ng/mL	ng/mL
Decimal Precision/Precision:	X.X	X.X
Reaction Direction:	Positive	Positive
Calculation Factor:	0	1
Math Model:	1	1
Cal. Time Limit:	336	336
Number of Calibrators:	5	5
#1	0.0	0.0
#2	250.0	250.0
#3	500.0	500.0
#4	750.0	750.0
#5	1000.0	1000.0
#6		
Primary Wavelength:	340 nm	340 nm
Secondary Wavelength:	650 nm	650 nm
Sample Volume:	15 µL	10 µL

REAGENTS:

Primary Inject (first)/First Inject		
Compartment/Component:	A	A
Volume/Dispense Volume:	200 µL	125 µL
Add Time/Inject Time:	NA ^a	NA ^a
Primary Inject (first)/Second Inject		
Compartment/Component:	None	None
Volume/Dispense Volume:		NA ^a
Add Time/Inject Time:	NA ^a	-180 ^{a&b}
Second Inject/Third Inject		
Compartment/Component:	B	B
Volume/Dispense Volume:	75 µL	75 µL
Add Time/Inject Time:	368 sec	276 sec

REAGENTS:

Blank		
Start Read:	237 sec	292 sec

Continued on next page

INSTRUMENT PARAMETERS:	SYNCHRON CX	UniCel DxC Synchron LX
End Read:	300 sec	308 sec
Reaction 1		
Start Read:	96 sec	336 sec
End Read:	144 sec	396 sec
Reaction 2		
Start Read:	NA ^a	NA ^a
End Read:	NA ^a	NA ^a
USABLE RESULT RANGE:		
Lower Limit:	0.00	0.00
Upper Limit:	1000.0	1000.0
ERROR DETECTION LIMITS:		
Reagent Blank/Blank		
ABS Low Limit:	-1.5 ^b	-1.5 ^b
ABS High Limit:	1.5 ^b	2.2 ^b
Rate Low Limit:	NA ^a	-1.5 ^b
Rate High Limit:	NA ^a	2.200 ^b
Mean Deviation:	NA ^a	2.200 ^b
Reaction/Reaction 1		
ABS Low Limit:	-1.5 ^b	-1.5 ^b
ABS High Limit:	1.5 ^b	2.2 ^b
Rate Low Limit:	NA ^a	-1.5 ^b
Rate High Limit:	NA ^a	2.200 ^b
Mean Deviation:	NA ^a	2.200 ^b
Reaction 2		
ABS Low Limit:	NA ^a	-1.5 ^b
ABS High Limit:	NA ^a	2.2 ^b
Rate Low Limit:	NA ^a	-1.5 ^b
Rate High Limit:	NA ^a	2.200 ^b
Mean Deviation:	NA ^a	2.200 ^b
SUBSTRATE DEPLETION		
Initial Rate:	99.999 ^b	99.999 ^b
Delta ABS:	1.5 ^b	2.2 ^b
MULTI POINT SPAN:		
(1-2)	0.004	0.009
(2-3)	0.024	0.046
(3-4)	0.031	0.037
(4-5)	0.016	0.013
(5-1)	0.082	0.106

^a NA=Not applicable

^b Denotes default value

Results and Data Interpretation

Refer to the package insert for information on data interpretation.

Typical Precision A properly operating Synchron system should exhibit precision values less than or comparable to the following:

Within-run Precision (Qualitative)

Rate mA/min	CX4			DxC600			LX20		
	Mean	SD	%CV	Mean	SD	%CV	Mean	SD	%CV
Negative Control (375 ng/mL)	83	0.76	0.91	115.6	1.50	1.29	118.3	1.77	1.49
Cutoff Calibrator (500 ng/mL)	102	0.83	0.81	146.2	1.95	1.33	149.6	1.91	1.28
Positive Control (625 ng/mL)	123	0.73	0.60	172.9	1.42	0.82	176.9	2.16	1.22

Total Run Precision (Qualitative)

Rate mA/min	CX4			DxC600			LX20		
	Mean	SD	%CV	Mean	SD	%CV	Mean	SD	%CV
Negative Control (375 ng/mL)	83	1.00	1.20	116	2.45	2.12	118	3.96	3.34
Cutoff Calibrator (500 ng/mL)	102	1.19	1.16	146	3.36	2.20	150	3.92	2.62
Positive Control (625 ng/mL)	123	1.08	0.89	173	3.08	1.78	177	4.25	2.40

Within-run Precision (Semi-quantitative)

Conc. ng/mL	CX4			DxC600			LX20		
	Mean	SD	%CV	Mean	SD	%CV	Mean	SD	%CV
Negative Control (375 ng/mL)	342.4	4.83	1.33	361.7	7.64	2.11	368.6	6.68	1.81
Cutoff Calibrator (500 ng/mL)	498.2	5.04	1.01	502.6	8.01	1.59	502.3	10.17	2.03
Positive Control (625 ng/mL)	633.5	6.67	1.05	634.8	13.07	2.06	629.7	11.49	1.82

Continued to next page

Total Precision (Semi-quantitative)

Conc. ng/mL	CX4			DxC600			LX20		
	Mean	SD	%CV	Mean	SD	%CV	Mean	SD	%CV
Negative Control (375 ng/mL)	342.4	8.86	2.45	361.7	9.24	2.55	368.6	13.24	3.59
Cutoff Calibrator (500 ng/mL)	498.2	7.09	1.42	502.6	9.80	1.95	502.3	10.61	2.11
Positive Control (625 ng/mL)	633.5	7.76	1.22	634.8	15.25	2.40	629.7	15.32	2.43

Continued on next page

**Cross
Reactivity**

Ecstasy compounds and ecstasy structurally related compounds were tested for cross reactivity. The table below presents representative results for Synchron instruments.

Ecstasy Compounds That Tested Positive

Cross-Reactant	Tested Conc. ng/mL
Cutoff Calibrator	
3,4-Methylenedioxyamphetamine (MDMA)	500
3,4-Methylenedioxyamphetamine (MDA)	900
3,4-Methylenedioxyethylamphetamine (MDEA)	400
N-Methylbenzodioxazolylbutananamine (MBDB)	2000
Benzodioxazolylbutananamine (BDB)	1500
p-Methoxyamphetamine (PMA)	4700
p-Methoxymethamphetamine (PMMA)	1700

Ecstasy Structurally Related Compounds That Tested Negative

Instruments	Synchron CX	UniCel DxC Synchron LX
Cross-Reactant Concentration	ng/mL	ng/mL
Cutoff Calibrator		
3,4-Methylenedioxyamphetamine (MDMA)	500	500
d-Amphetamine	600,000	600,000
l-Amphetamine	10,000	30,000
d,l-Amphetamine	40,000	75,000
l-Ephedrine	800,000	800,000
d-Methamphetamine	200,000	300,000
l-Methamphetamine	10,000	10,000
d,l-Methamphetamine	20,000	30,000
Phentermine	50,000	100,000
d,l-Phenylpropanolamine	800,000	800,000
d-Pseudoephedrine	1,000,000	1,000,000

Continued on next page

Method Comparison

Refer to the Accuracy Section of the Product Insert for GC/MS^c Comparison

A total of 138 clinical samples were tested in the **Qualitative and Semi-quantitative** modes on the Synchron and compared to the results by GC/MS. The data were analyzed and presented in the table below.

Qualitative

CX4				DxC600				LX20			
		+	-			+	-			+	-
GC/MS	+	50	0	GC/MS	+	50	0	GC/MS	+	50	0
	-	7 ^c	81		-	7 ^c	81		-	7 ^c	81
Positive Sample Agreement			100%	Positive Sample Agreement:			100%	Positive Sample Agreement:			100%
Negative Sample Agreement			92%	Negative Sample Agreement			92%	Negative Sample Agreement			92%
Total Sample Agreement			94.9%	Total Sample Agreement			94.9%	Total Sample Agreement			94.9%

Semi-quantitative:

CX4				DxC600				LX20			
		+	-			+	-			+	-
GC/MS	+	50	0	GC/MS	+	50	0	GC/MS	+	50	0
	-	7 ^c	81		-	7 ^c	81		-	7 ^c	81
Positive Sample Agreement			100%	Positive Sample Agreement:			100%	Positive Sample Agreement:			100%
Negative Sample Agreement			92%	Negative Sample Agreement			92%	Negative Sample Agreement			92%
Total Sample Agreement			94.9%	Total Sample Agreement			94.9%	Total Sample Agreement			94.9%

^c Of the 138 total clinical samples confirmed by GC/MS (Gas Chromatography / Mass Spectrometry) seven were between 500 ng/ml and 634 ng/ml (+25% of cutoff concentration). Due to limitations in sample volume, not all Ecstasy derivatives and/or metabolites were tested for.

ADDITIONAL INFORMATION

This guideline has undergone limited technical evaluation and is intended to provide guidance only for the use of this reagent on the UniCel[®] Dx_C, SYNCHRON LX[®] and CX[®] Clinical Systems. The performance characteristics have not been established. You should perform additional testing before reporting diagnostic results.

Information on sample preparation, expected values, quality control, as well as warnings and precautions related to the use of this reagent may be obtained from the package insert.

Instrument operating instructions are contained in the SYNCHRON LX[®] Operations Manual, SYNCHRON CX[®] Operating Instructions, or UniCel[®] Dx_C Systems Instructions For Use (IFU) Manual.

Since Beckman Coulter does not manufacture the reagent or perform quality control or other tests on individual lots, Beckman Coulter cannot be responsible for the quality of the data obtained which is caused by performance of the reagent, any variation between lots of reagent, or protocol changes by the manufacturer.

SHIPPING DAMAGE

If damaged product is received, notify your Beckman Coulter Clinical Support Center.