

DRI[®] TRICYCLICS SERUM TOX APPLICATION
Beckman Coulter AU400[®], AU480[®], AU640[®], AU680[®],
AU2700[®], AU5400[®], AU5800[®]



Catalog No. 1128

The DRI Tricyclics Serum Tox Assay is intended for the qualitative and semiquantitative determination of tricyclic antidepressants in human serum or plasma or 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	Thermo Fisher Scientific Catalog Number
DRI Tricyclics Serum Tox Kit	R1 25 mL R2 8 mL	1128
DRI Serum Tox Negative Calibrator	10 mL	0962
DRI Serum Tox Calibrator 1	5 mL	0963
DRI Serum Tox Calibrator 2	5 mL	0965
DRI Serum Tox Calibrator 3	5 mL	0967
DRI Serum Tox Calibrator 4	5 mL	0976
MAS TOX Control 1 – 3	6 x 5 mL	10011608

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

Microgenics GmbH, Spitalhofstrasse 94, 94032 Passau, Germany

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.

**Results and
Data
Interpretation**

Results for samples will be printed in ng/mL.

DRI Tricyclics Serum Tox Qualitative Application

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

Specific Test Parameters	
General	LIH ISE Range
Test Name:	TCACQ < >
Type:	Serum <
Operation:	Yes <
Sample: Volume	8.0 μL
Dilution	0 μL
Pre-Dilution Rate:	1
Reagents: R1 Volume	180 μL
Dilution	0 μL
Min OD	
Max OD	
R2 Volume	60 μL
Dilution	0 μL
Reagent OD limit:	L -2.00 H 2.50
Wavelength: Pri.	340 <
Sec.	520 <
First L	-2.00
First H	2.50
Method:	FIXED1* <
Last L	-2.00
Last H	2.50
Reaction slope:	+ <
Dynamic Range:	L # H #
Measuring Point 1: First	14
Last	20
Measuring Point 2: First	
Last	
Correlation Factor:	A 1 B 0
Linearity	%
On-board stability period:	#
No-Lag-Time:	No <

Specific Test Parameters	
General	LIH ISE Range
Test Name:	TCACQ < >
Type:	Serum <
Value/Flag:	# <
Level L:	#
Level H:	†
Normal Ranges:	Age L Age H
	Sex Year Month Year Month L H
<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: % Decimal places: #

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

User defined
 * Can also be run as RATE1

DRI Tricyclics Serum Tox Semiquantitative Application

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

Specific Test Parameters	
General	LIH ISE Range
Test Name:	TCASQ < > Type: Serum Operation: Yes
Sample:	Volume: 8.0 μL Dilution: 0 μL Pre-Dilution Rate: 1
Reagents:	R1 Volume: 180 μL Dilution: 0 μL Min OD Max OD
	R2 Volume: 60 μL Dilution: 0 μL L: -2.00 H: 2.50
Wavelength:	Pri.: 340 Sec.: 520 Reagent OD limit: First L: -2.00 First H: 2.50
Method:	FIXED1* Last L: -2.00 Last H: 2.50
Reaction slope:	+
Measuring Point 1:	First: 14 Last: 20 L: 0 H: 1000
Measuring Point 2:	First: Last: Correlation Factor: A: 1.0000 B: 0.0000
Linearity:	% On-board stability period: #
No Lag Time:	No

Specific Test Parameters	
General	LIH ISE Range
Test Name:	TCASQ < > Type: Serum
Value/Flag:	# Level L: # Level H: #
Normal Ranges:	Age L: Age H: L: H:
	Sex Year Month Year Month L H
<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:	TCASQ < > Type: Serum
Calibration Type:	5AB Formula: EIA TYPE 1 Counts: 2 Process: CONC
Point 1:	Cal. No.: # OD: CONC: 0 Factor/OD-L: -2.000 Factor/OD-H: 2.5000
Point 2:	# # # -2.000 2.5000
Point 3:	# # # -2.000 2.5000
Point 4:	# # # -2.000 2.5000
Point 5:	# # # -2.000 2.5000
Point 6:	# # # # #
Point 7:	# # # # #
1-Point Cal. Point:	# <input type="checkbox"/> With CONC-0 Slope Check + Advanced Calibration: #
MB Type Factor:	Calibration Stability Period: #

* Can also be run as RATE1

User defined

DRI Tricyclics Serum Tox Qualitative Application

Beckman Coulter System Parameters, AU480/AU680

Specific Test Parameters

General LIH ISE Range

Test Name: TCACQ < > Type: Serum Operation: Yes

Sample Volume: 7.2 μL Dilution: 0 μL OD Limit: Min. OD: -2.00 Max. OD: 3.00

Pre-Dilution Rate: 1

Reagents Volume: R1(R1-1) 162 μL Dilution: 0 μL Reagent OD limit: First Low: -2.00 High: 3.00 Last Low: -2.00 High: 3.00

R2 (R2-1) 54 μL Dilution: 0 μL Dynamic Range Low: # High: #

Correlation Factor A: 1 B: 0

Factor for Maker A: 1 B: 0

Wavelength: Pri. 340 nm Sec. 520 nm

Method: FIXED1*

Reaction slope: +

Onboard Stability: 32 Days 0 Hour

Measuring Point 1: First 14 Last 20 LIH Influence Check: #

Measuring Point 2: First Last

Linearity: %

No Lag Time: No

Lipemia: Icterus: Hemolysis:

Specific Test Parameters

General ISE Range

Test Name: TCACQ < > Type: Serum

Value/Flag: # Level L: # Level H: †

Specific Ranges:

	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: #

Panic Value: Low: # High: #

Calibration Specific

General ISE

Test Name: TCACQ < > Type: Urine Use Serum Cal.

Calibration Type: AA Formula: Y=AX+B Counts: 2

Slope Check: +

Calibrator †	OD	Conc	Factor Range		Allowable Range Check
			Low	High	
Point 1: #		0	-99999	999999	<input type="checkbox"/> Reagent Blank
Point 2: #		100	-99999	999999	<input type="checkbox"/> Calibration
Point 3:					
Point 4:					
Point 5:					
Point 6:					
Point 7:					
Point 8:					
Point 9:					
Point 10:					

Advanced Calibration: Operation Interval (RB/ACAL):

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

Calibrator	OD	Conc	OD Range		Stability
			Low	High	
Point 1:					# Day # Hour
Point 2:					18 Day 0 Hour

Reagent Blanks Calibration: # Day # Hour

MB Type Factor: 1-Point Calibration Point: With CONC-0:

User defined.

* Can also be run as RATE1

DRI Tricyclics Serum Tox Semiquantitative Application

Beckman Coulter System Parameters, AU480/AU680

Specific Test Parameters														
General		LIH	ISE	Range										
Test Name:	TCASQ	<	>	Type:	Serum	Operation:	Yes							
Sample Volume	7.2	μL	Dilution	0	μL	OD Limit								
Pre-Dilution Rate	1					Min. OD	-2.000	Max. OD	3.00					
Reagents Volume:	R1(R1-1)	162	μL	Dilution	0	μL	Reagent OD limit:							
							First Low	-2.000	High	3.000				
							Last Low	-2.000	High	3.000				
	R2 (R2-1)	54	μL	Dilution	0	μL	Dynamic Range Low	0	High	1000				
Wavelength:	Pri.	340	nm	Sec.	520	nm	Correlation Factor A	1	B	0				
Method:	FIXED1*													
Reaction slope:	+													
Measuring Point 1:	First	14	Last	20	Onboard Stability						32	Days	0	Hour
Measuring Point 2:	First		Last		LIH Influence Check						#			
Linearity:											Lipemia			
No Lag Time:	No										Icterus			
											Hemolysis			

Specific Test Parameters										
General		ISE	Range							
Test Name:	TCASQ	<	>	Type:	Serum					
Value/Flag:	#	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	#						

Calibration Specific										
General		ISE								
Test Name:	TCASQ	<	>	Type:	Serum	<input type="checkbox"/> Use Serum Cal.				
Calibration Type:	5AB	Formula:	POLYGONAL	Counts:	2	Slope Check	+			
<Calibrator Parameters>										
Calibrator ↑	OD	Conc	Factor Range		Allowable Range Check					
			Low	High	<input type="checkbox"/> Reagent Blank					
Point 1:	#	0	-2.0000	3.0000	<input type="checkbox"/> Calibration					
Point 2:	#	150	-2.0000	3.0000	Advanced Calibration					
Point 3:	#	300	-2.0000	3.0000	Operation					
Point 4:	#	500	-2.0000	3.0000	Interval (RB/ACAL)					
Point 5:	#	1000	-2.0000	3.0000						
Point 6:										
Point 7:										
Point 8:										
Point 9:										
Point 10:										
<Point Cal. For Master Curve>										
Calibrator	OD	Conc	OD Range		Stability					
			Low	High	Reagent Blanks					
Point 1:					#					
Point 2:					Day					
					#					
					Hour					
MB Type Factor: # 1-Point Calibration Point # With CONC-0										

User Defined
 * Can also be run as RATE1

DRI Tricyclics Serum Tox Qualitative Application

Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters									
General	LIH	ISE	HbA1c		Calculated Test	Range					
Test Name:		TCACQ	<	>	Type:	Serum	Operation	Yes			
Sample Volume	7.2	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)	162	μL	Dilution	0	μL	Reagent OD Limit				
	R1-2		μL	Dilution		μL	1 st .	Low	-2.00	High	3.00
	R2(R2-1)	54	μL	Dilution	0	μL	Last	Low	-2.00	High	3.00
Common Rgt. Type	None		Name			Correlation Factor A	1				
Wavelength	Pri	340	▽nm	Sec.	520	▽nm	Factor for Maker A	1			
Method	FIXED1*						High	#			
Reaction Slope	+						B	0			
Measuring Point1 1 st	14		Last	20		Onboard Stability Period	32	Day	0	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 Test	Range		
Test Name:		TCACQ	<	>	Type:	Serum	▽	
Value/Flag:	#		Level		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	%	Decimal Places	#

Parameters		Calibration Parameters						
Calibrators	Calibration Specific							
General	ISE							
Test Name:		TCACQ	<	>	Type:	Serum	Cuvette .	▽
								<input type="checkbox"/> Use Serum Cal.
Calibration Type:	AA	Formula:	Y=AX+B		Counts:	2		
<Calibrator Parameters>		Range		Slope Check		+		
Calibrator	OD	Conc	Low	High	Allowance Range Check			
Point 1:	#	0	-99999	999999	<input type="checkbox"/> Reagent Blank			
Point 2:	#	100	-99999	999999	<input type="checkbox"/> Calibration			
Point 3:					Advanced Calibration			
Point 4:					Operation			
Point 5:					No		▽	
Point 6:					Interval (RB/ACAL)			
Point 7:								
Point 8:								
Point 9:								
Point 10:								
<Point Cal. For	No. of Correction Points		Use Master Curve		<input type="checkbox"/> Lot Calibration			
Master Curve>		OD Range		Stability				
Calibrator	OD	Conc	Low	High	Reagent Blank	#	Day	#
Point-1					Calibration	18	Day	0
Point-2								
MB Type Factor:	1-Point Calibration Point		None		<input type="checkbox"/> with Conc-0			

User defined

* Can also be run as RATE1

DRI Tricyclics Serum Tox Semiquantitative Application

Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: TCASQ				Type: Serum	Operation: Yes
Sample Volume	7.2	μL	Dilution	0	μL
Pre-Dilution Rate	1		Diluent Bottle	#	
Rgt. Volume	R1(R1-1)	162	μL	Dilution	0
	R1-2		μL	Dilution	
	R2(R2-1)	54	μL	Dilution	0
Common Rgt. Type	None		Name	None	
Wavelength	Pri	340	nm	Sec.	520
Method	FIXED1*		Dynamic Range	Low	0
Reaction Slope	+		Correlation Factor	A	1
Measuring Point1 1 st	14		Factor for Maker	A	1
Measuring Point2 1 st			High	1000	
Linearity Limit			Low	-2.00	
Lag Time Check	No		High	3.00	
			Onboard Stability Period	32 Day	
			LIH Influence Check	#	
			Lipemia		
			Icterus		
			Hemolysis		

Parameters		Specific Test Parameters			
General	LIH	ISE	HbA1c	Calculated Test	Range
Test Name: TCASQ				Type: Serum	
Value/Flag: #					
Specific Ranges:		From	Level 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. Standard demographics			#	#
	8. Not within expected values			#	#
Panic Value	Low	#	High	#	Unit: ng/mL
					Decimal Places: #

Parameters		Calibration Parameters			
Calibrators	Calibration Specific				
General	ISE				
Test Name: TCASQ				Type: Serum	Cuvette: #
		<input type="checkbox"/> Use Serum Cal.			
Calibration Type:	5AB	Formula:	POLYGONAL		Counts: 2
<Calibrator Parameters>		Calibrator	OD	Conc	Range
Point 1:	#			0	Low: -2.00 High: 3.00
Point 2:	#			150	Low: -2.00 High: 3.00
Point 3:	#			300	Low: -2.00 High: 3.00
Point 4:	#			500	Low: -2.00 High: 3.00
Point 5:	#			1000	Low: -2.00 High: 3.00
Point 6:					
Point 7:					
Point 8:					
Point 9:					
Point 10:					
<Point Cal. For Master Curve>	No. of Correction Points		Use Master Curve		<input type="checkbox"/> Lot Calibration
Calibrator	OD	Conc	Low	High	Stability
Point-1					Reagent Blank # Day # Hour
Point-2					Calibration 18 Day 0 Hour
MB Type Factor:	1-Point Calibration Point		None		<input type="checkbox"/> with Conc-0

User defined
 * Can also be run as RATE1

Precision

Tests for within-run and total precision, evaluated with packaged reagents, controls and calibrators, yielded the following results (N=80/level):

Controls	Control 1	Control 2	Control 3	Cutoff
Qualitative AU480				
Mean (mA/min)	796	870	923	837
Within-Run SD (mA/min)	3.0	3.0	2.6	2.8
Within-Run CV (%)	0.4%	0.3%	0.3%	0.3%
Total SD (mA/min)	3.8	3.6	3.3	4.2
Total CV (%)	0.5%	0.4%	0.4%	0.5%
Qualitative AU640*				
Mean (mA/min)	454	590	N/A	536
Within-Run SD (mA/min)	3.1	3.6	N/A	4.4
Within-Run CV (%)	0.7%	0.6%	N/A	0.8%
Total SD (mA/min)	4.1	4.8	N/A	5.2
Total CV (%)	0.9%	0.8%	N/A	1.0%
Qualitative AU680				
Mean (mA/min)	786	862	913	833
Within-Run SD (mA/min)	5.7	5.7	6.1	6.2
Within-Run CV (%)	0.7%	0.7%	0.7%	0.7%
Total SD (mA/min)	31.6	27.5	31.8	24.8
Total CV (%)	4.0%	3.2%	3.5%	3.0%
Qualitative AU5800				
Mean (mA/min)	771	844	897	811
Within-Run SD (mA/min)	11.4	12.3	12.0	10.9
Within-Run CV (%)	1.5%	1.5%	1.3%	1.3%
Total SD (mA/min)	11.5	12.5	12.3	11.2
Total CV (%)	1.5%	1.5%	1.4%	1.4%

Controls	Control 1	Control 2	Control 3	Cutoff
Semiquantitative AU480				
Mean (ng/mL)	225	439	786	300
Within-Run SD (ng/mL)	5.2	12.9	19.0	7.8
Within-Run CV (%)	2.3%	2.9%	2.4%	2.6%
Total SD (ng/mL)	6.0	15.7	25.7	9.7
Total CV (%)	2.7%	3.6%	3.3%	3.2%
Semiquantitative AU640*				
Mean (ng/mL)	158	548	N/A	317
Within-Run SD (ng/mL)	6.0	31.0	N/A	14.2
Within-Run CV (%)	3.8%	5.7%	N/A	4.5%
Total SD (ng/mL)	7.8	39.4	N/A	16.9
Total CV (%)	4.9%	7.2%	N/A	5.3%
Semiquantitative AU680				
Mean (ng/mL)	218	434	808	302
Within-Run SD (ng/mL)	9.1	29.3	50.6	17.0
Within-Run CV (%)	4.2%	6.7%	6.3%	5.6%
Total SD (ng/mL)	16.5	41.4	63.2	19.8
Total CV (%)	7.5%	9.5%	7.8%	6.6%
Semiquantitative AU5800				
Mean (ng/mL)	224	436	777	300
Within-Run SD (ng/mL)	11.3	23.6	40.2	12.6
Within-Run CV (%)	5.1%	5.4%	5.2%	4.2%
Total SD (ng/mL)	12.1	26.3	45.7	14.8
Total CV (%)	5.4%	6.0%	5.9%	4.9%

AU640* N=60/level

Accuracy and Correlation AU480, AU680, and AU5800

Patient samples were assayed using the DRI Tricyclics Serum Tox Assay on the Beckman Coulter AU480, AU680, and AU5800, and tested against the reference analyzer, the Hitachi 717.

Qualitative Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	90.2%	100.0%	96.4%
AU680	97.6%	97.1%	97.3%
AU5800	95.8%	98.9%	97.9%

DRI Tricyclics Serum Tox Assay Qualitative Method Comparison											
		Hitachi 717				Hitachi 717				Hitachi 717	
		+	-			+	-			+	-
AU480	+	46	0	AU680	+	41	2	AU5800	+	46	1
	-	5	89		-	1	66		-	2	91

Semiquantitative Method Comparison			
	Positive Agreement (%)	Negative Agreement (%)	Total Agreement (%)
AU480	97.8%	96.8%	97.1%
AU680	93.5%	100.0%	97.3%
AU5800	95.7%	100.0%	98.2%

DRI Tricyclics Serum Tox Assay Semiquantitative Method Comparison											
		Hitachi 717				Hitachi 717				Hitachi 717	
		+	-			+	-			+	-
AU480	+	44	3	AU680	+	43	0	AU5800	+	44	0
	-	1	92		-	3	64		-	2	64

AU640 Accuracy and Correlation

The AU640 data is presented as being representative of the performance of this instrument family.

One hundred twenty seven (127) serum samples were assayed in qualitative mode using the DRI Tricyclics Serum Tox Assay on the Beckman AU640 and the Hitachi 717 analyzers. A Sensitivity of 100% (89 of 89 positive samples) and a Specificity of 98% (50 of 51 negative samples) were observed between the two analyzers.

A linear regression analysis of the semi-quantitative data yielded the following results:
 Beckman AU640 = 1.04 (Hitachi 717) – 6.7, with a correlation coefficient of 0.950

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

End