

CEDIA® TOBRAMYCIN II APPLICATION



Beckman Coulter AU400® , AU480® , AU640® , AU680® , AU2700® , AU5400® , AU5800®

Catalog No. 100018

Homogeneous Enzyme Immunoassay for the Quantitative Determination of Tobramycin Levels in Serum and Plasma

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 No
CEDIA Tobramycin II Assay Reagents	13 mL 11 mL	100018
CEDIA TDM Antibiotic Multi-Cal	2x5 mL 2x7.5 mL	100017
MAS PAR TDM Level 1	6x5 mL	PTD1-1001
MAS PAR TDM Level 2	6x5 mL	PTD2-2002
MAS PAR TDM Level 3	6x5 mL	PTD3-3003

To place an order or for technical service contact:

USA	In Europe
(800) 232-3342 Fax 510-979-5002	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.

Analyzer Procedure

Refer to the operator's manuals for information on analyzer operation.

Dispense adequate amounts of EA Reagent (first reagent) and ED Reagent (second reagent) into appropriate containers place reagents onboard the analyzer prior starting analysis. **Ensure that the reagents have equilibrated to the temperature of the analyzer reagent compartment before starting analysis**

Note: Under Specific Test Parameters/General Tab, Linearity % should be left blank, as reflected in the following pages. **Do Not Enter 0.**

Calibration

Refer to the package insert for information on calibration.

Results and Data Interpretation

For further information, contact Microgenics Technical Support or your local Microgenics representative.

CEDIA TOBRAMYCIN Application
Beckman Coulter System Parameters, AU400/AU640/AU2700/AU5400

Specific Test Parameters											
General		LIH	ISE	Range							
Test Name:		TOBRA	<	>	Type:	Serum	Operation:	Yes			
Sample Volume		6.5	μL	Dilution	0	μL	OD Limit				
Pre-Dilution Rate		1					Min. OD	-2.00	Max. OD	2.50	
Reagents Volume:	R1(R1-1)	82	μL	Dilution	0	μL	Reagent OD limit:				
							First Low	-2.00	High	2.50	
							Last Low	-2.00	High	2.50	
	R2 (R2-1)	62	μL	Dilution	0	μL	Dynamic Range Low	*	High	*	
							Correlation Factor A	1	B	0	
Wavelength:	Pri.	570	nm	Sec.	660	nm	Factor for Maker A		B		
Method:		RATE1									
Reaction slope:		+					Onboard Stability	#	Days	0	Hour
Measuring Point 1:	First	24		Last	27		LIH Influence Check	#			
Measuring Point 2:	First			Last			Lipemia				
Linearity:			%				Icterus				
No Lag Time:		No					Hemolysis				

Specific Test Parameters											
General		ISE	Range								
Test Name:		TOBRA	<	>	Type:						
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	*	Decimal Places					#				

CEDIA TOBRAMYCIN Assay
Beckman Coulter System Parameters, AU400/AU640/AU2700/AU5400, Continued

Calibration Specific										
General		ISE								
Test Name:		TOBRA		<	>	Type				<input type="checkbox"/> Use Serum Cal.
Calibration Type:		2AB		Formula:		POLYGONAL		Counts:	2	
<Calibrator Parameters>										
	Calibrator ↑	OD	Conc	Factor Range				Slope Check	+	
				Low	High			Allowable Range Check		
Point 1:	#		**	-2.00	3.00			<input type="checkbox"/> Reagent Blank		
Point 2:			**	-2.00	3.00			<input type="checkbox"/> Calibration		
Point 3:								Advanced Calibration		
Point 4:								Operation		
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	Day	Hour
Point 1:								Calibration	Day	Hour
Point 2:										
MB Type Factor:				1-Point Calibration Point				<input type="checkbox"/> With CONC-0		

User Defined
 ** Lot specific calibrator value.

CEDIA TOBRAMYCIN Assay
Beckman Coulter System Parameters, AU480 & AU680

Specific Test Parameters															
General		LIH	ISE	Range											
Test Name:		TOBRA	<	>	Type:	Serum	Operation:	Yes							
Sample Volume	6.5	μL	Dilution	0	μL	OD Limit									
Pre-Dilution Rate	1					Min. OD	-2.00	Max. OD	3.00						
Reagents Volume:	R1(R1-1)	82	μL	Dilution	0	μL	Reagent OD limit:								
						First Low	-2.00	High	3.00						
						Last Low	-2.00	High	3.00						
	R2 (R2-1)	62	μL	Dilution	0	μL	Dynamic Range Low	0.07	High	11.8					
						Correlation Factor A	1	B	0						
Wavelength:	Pri.	570	nm	Sec.	660	nm	Factor for Maker A		B						
Method:	FIXED1* ▾														
Reaction slope:	+		▾								Onboard Stability	30	Days	0	Hour
Measuring Point 1:	First	24	Last	27	LIH Influence Check								#	▾	
Measuring Point 2:	First		Last		Lipemia									▾	
Linearity:											Icterus		▾		
No Lag Time:	No		▾								Hemolysis		▾		

Specific Test Parameters											
General		ISE	Range								
Test Name:		TOBRA	<	>	Type:						
Value/Flag:	#	▾	Level L:	#	Level H:	#					
Specific Ranges:										Panic Value	
	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	#

* Can also be run as RATE1

User Defined Values

CEDIA TOBRAMYCIN
Beckman Coulter System Parameters, AU480 & AU680, Continued

Calibration Specific										
General		ISE								
Test Name:		TOBRA ▾		<	>	Type		▾		<input type="checkbox"/> Use Serum Cal.
Calibration Type:		2AB ▾		Formula:		POLYGONAL ▾		Counts:		#
<Calibrator Parameters>										
Calibrator ↑		OD	Conc	Factor Range		Slope Check		+		▾
				Low	High	Allowable Range Check				
Point 1:	# ▾		**	-2.00	3.00	<input type="checkbox"/> Reagent Blank				
Point 2:	# ▾		**	-2.00	3.00	<input type="checkbox"/> Calibration				
Point 3:	▾					Advanced Calibration				
Point 4:	▾					Operation		▾		
Point 5:	▾					Interval (RB/ACAL)		▾		
Point 6:	▾									
Point 7:	▾									
Point 8:	▾									
Point 9:	▾									
Point10:	▾									
<Point Cal. For Master Curve>		No. of Correction Points		▾		Use Master Curve		▾		<input type="checkbox"/> Lot Calibration
Calibrator		OD	Conc	OD Range		Stability				
				Low	High	Reagent Blanks Calibration				
Point 1:						3 Day				Hour
Point 2:						0 Day				Hour
MB Type Factor:		1-Point Calibration Point		▾		<input type="checkbox"/> With CONC-0				

User Defined
 ** Lot specific calibrator value.

CEDIA TOBRAMYCIN Assay
Beckman Coulter System Parameters, AU5800

Parameters		Specific Test Parameters									
General	LIH	ISE	HbA1c		Calculated Test	Range					
Test Name:		TOBRA	<	>	Type:	Urine	Operation	Yes			
Sample Volume	6.5	μL	Dilution	0	μL	OD Limit					
Pre-Dilution Rate	1	∇	Diluent Bottle	#	∇	Min.OD	-2.000	Max.OD	3.000		
Rgt. Volume	R1(R1-1)	82	μL	Dilution	0	μL	Reagent OD Limit				
	R1-2		μL	Dilution		μL	1 st .	Low	-2.000	High	3.000
							Last	Low	-2.000	High	3.000
	R2(R2-1)	62	μL	Dilution	0	μL					
Common Rgt. Type	None		Name	None		Dynamic Range Low	0.07	High	11.8		
Wavelength	Pri	570	∇nm	Sec.	660	∇nm	Correlation Factor A	1	B	0	
Method	FIXED1*					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						Hemolysis		∇			

Parameters		Specific Test Parameters							
General	LIH	ISE	HbA1c		Calculated Test	Range			
Test Name:		TOBRA	<	>	Type:	Urine			
Value/Flag:		#							
Specific Ranges:		From		Level To		Low	#	High	#
	Sex	Year	Month	Year	Month	Low		High	
o 1.	#	#	#	#	#	#	#	#	#
o 2.	#	#	#	#	#	#	#	#	#
o 3.	#	#	#	#	#	#	#	#	#
o 4.	#	#	#	#	#	#	#	#	#
o 5.	#	#	#	#	#	#	#	#	#
o 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:		TOBRA	<	>	Type	Urine	Cuvette .	
		<input type="checkbox"/> Use Serum Cal.						
Calibration Type:		2AB	Formula:	Polygonal	Counts:	#		
<Calibrator Parameters>		Range		Slope Check		+		
Calibrator	OD	Conc	Low	High				
Point 1:	#	**	-2.000	3.000	Allowance Range Check			
Point 2:	#	**	-2.000	3.000	<input type="checkbox"/> Reagent Blank			
Point 3:					<input type="checkbox"/> Calibration			
Point 4:					Advanced Calibration			
Point 5:					Operation			
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	Hour	
Point-1					3	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.
 ** Lot specific calibrator value.

Precision

Three control levels were tested in replicates of 2, twice per day for 20 days, total N = 80. The results are presented in the following tables:

*N=60

	C1	C2	C3
Quantitative AU400*			
Mean (ng/mL)	1.8	4.7	7.8
Within-Run SD (ng/mL)	0.11	0.16	0.23
Within-Run CV (%)	6.3	3.5	3.0
Total SD (ng/mL)	0.16	0.19	0.26
Total CV (%)	9.1	4.0	3.4
Quantitative AU480			
Mean (ng/mL)	2.0	4.6	7.9
Within-Run SD (ng/mL)	0.04	0.06	0.19
Within-Run CV (%)	2.1	1.3	2.4
Total SD (ng/mL)	0.07	0.09	0.34
Total CV (%)	3.3	2.0	4.2
Quantitative AU640*			
Mean (ng/mL)	1.7	4.7	7.8
Within-Run SD (ng/mL)	0.08	0.13	0.17
Within-Run CV (%)	4.8	2.9	2.2
Total SD (ng/mL)	0.10	0.18	0.21
Total CV (%)	5.7	4.0	2.7
Quantitative AU680			
Mean (ng/mL)	2.0	4.6	8.0
Within-Run SD (ng/mL)	0.07	0.10	0.13
Within-Run CV (%)	3.6	2.1	1.6
Total SD (ng/mL)	0.08	0.13	0.18
Total CV (%)	3.9	2.8	2.2
Quantitative AU5800			
Mean (ng/mL)	2.0	4.6	8.0
Within-Run SD (ng/mL)	0.09	0.08	0.16
Within-Run CV (%)	4.2	1.8	2.0
Total SD (ng/mL)	0.12	0.16	0.24
Total CV (%)	5.9	3.5	3.0

**Assay
Accuracy and
Correlation
Beckman
AU400**

One hundred twenty-four (124) samples were assayed with the Tobramycin II Assay on both the Beckman AU400 and Hitachi 911 analyzers with the following results:
 $y = 0.98x + 0.06$; $r = 0.999$

**Assay
Accuracy and
Correlation
Beckman
AU480**

One hundred and eleven (111) samples were assayed with the Tobramycin II Assay on both the Beckman AU480 and Hitachi 911 analyzers with the following results:
 $y = 1.02x + 0.22$; $r = 0.998$

**Assay
Accuracy and
Correlation
Beckman
AU640**

One hundred twenty-four (124) samples were assayed with the Tobramycin II Assay on both the Beckman AU640 and Hitachi 911 analyzers with the following results:
 $y = 0.98x + 0.14$; $r = 0.997$

**Assay
Accuracy and
Correlation
Beckman
AU680**

One hundred and eleven (111) samples were assayed with the Tobramycin II Assay on both the Beckman AU680 and Hitachi 911 analyzers with the following results:
 $y = 1.00x + 0.24$; $r = 0.998$

**Assay
Accuracy and
Correlation
Beckman
AU5800**

One hundred and fifteen (115) samples were assayed with the Tobramycin II Assay on both the Beckman AU5800 and Hitachi 911 analyzers with the following results:
 $y = 1.05x + 0.22$; $r = 0.998$

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

End