

QMS™ TOBRAMYCIN APPLICATION
BECKMAN COULTER AU480® , AU680® , AU5800®

Catalog No. 10017109

This Application is Intended for the Quantitative Determination of Tobramycin in Human Serum



For In Vitro Diagnostic Use Only
Rx 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. For package inserts, visit www.thermofisher.com and enter the assay name in the *Search* field.

Ordering Information

For ordering information, visit www.thermofisher.com and enter the assay name in the *Search* field. Not all intended uses and applications mentioned here are available in every country. Please contact your local sales representative or distributor for more information.

For Technical Service, contact:

USA	In Europe
<p>Tel: (800) 232-3342 Fax: (510) 979-5420</p>	<p>Tel: +49 (0)851-88 6890 Fax: +49 (0)851-88 68910</p>



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 the package insert for information on reagent storage.

**Procedure for
Analyzer**

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

**Results and
Data
Interpretation**

Results for samples will be printed in $\mu\text{g/mL}$.

QMS Tobramycin Assay - Quantitative

Beckman Coulter System Parameters, AU480/AU680

Parameters		Specific Test Parameters									
General	LIH	ISE	HbA1c		Calculated Test	Range					
Test Name:		<input type="text" value="QTOBR"/>	<input type="text" value="<"/>	<input type="text" value=">"/>	Type:	<input type="text" value="#"/>	Operation	<input type="text" value="Yes"/>			
Sample Volume	<input type="text" value="6"/>	μL	Dilution	<input type="text" value="0"/>	μL	OD Limit					
Pre-Dilution Rate	<input type="text" value="1"/>	▽	Diluent Bottle	<input type="text"/>	▽	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="100"/>	μL	Dilution	<input type="text" value="0"/>	μL	Reagent OD Limit				
	R1-2	<input type="text"/>	μL	Dilution	<input type="text"/>	μL	1 st .	Low	<input type="text" value="-2.00"/>	High	<input type="text" value="3.00"/>
	R2(R2-1)	<input type="text" value="83"/>	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"/>		Dynamic Range Low	<input type="text" value="0.4"/>	High	<input type="text" value="10"/>		
Wavelength	Pri	<input type="text" value="700"/>	nm	Sec.	<input type="text"/>	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="+"/>	▽				Onboard Stability Period	<input type="text" value="#"/>	Day	<input type="text" value="#"/>	Hour	<input type="text" value="#"/>
Measuring Point1	1 st	<input type="text" value="14"/>		Last	<input type="text" value="19"/>		LIH Influence Check	<input type="text" value="#"/>	▽		
Measuring Point2	1 st	<input type="text"/>		Last	<input type="text"/>		Lipemia	<input type="text" value="++++"/>	▽		
Linearity Limit	<input type="text"/>	%					Icterus	<input type="text" value="++++"/>	▽		
Lag Time Check	<input type="text" value="No"/>	▽					Hemolysis	<input type="text" value="++++"/>	▽		

Parameters		Specific Test Parameters						
General	LIH	ISE	HbA1c		Calculated Test	Range		
Test Name:		<input type="text" value="#"/>	<input type="text" value="<"/>	<input type="text" value=">"/>	Type:	<input type="text" value="#"/>		
Value/Flag:		<input type="text" value="#"/>						
			Level		Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>
Specific Ranges:		From	To					
	Sex	Year	Month	Year	Month	Low	High	
<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="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="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="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="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="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="#"/>	<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="μg/mL"/>	Decimal Places	<input type="text" value="#"/>

User defined.

QMS Tobramycin Assay - Quantitative

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

Parameters		Calibration Parameters			
Calibrators		Calibration Specific			
General		ISE			
Test Name: QTOBR ▾		< > Type # ▾		Cuvette . ▾	
<input type="checkbox"/> Use Serum Cal.					
Calibration Type: 6AB ▾		Formula: EIA Type1 ▾		Counts: 2 ▾	
<Calibrator Parameters>					
	Calibrator	OD	Conc	Range	
Point 1:	# ▾		0†	Low	High
Point 2:	# ▾		0.5†	-2.00	3.00
Point 3:	# ▾		1.5†	-2.00	3.00
Point 4:	# ▾		3.0†	-2.00	3.00
Point 5:	# ▾		6.0†	-2.00	3.00
Point 6:	# ▾		10.0†	-2.00	3.00
Point 7:	# ▾				
Point 8:	# ▾				
Point 9:	# ▾				
Point 10:	# ▾				
				Slope Check - ▾	
				Allowance Range Check	
				<input type="checkbox"/> Reagent Blank	
				<input type="checkbox"/> Calibration	
				Advanced Calibration Operation ▾	
				Interval (RB/ACAL) ▾	
				<input type="checkbox"/> Lot Calibration	
<Point Cal. For Master Curve>		No. of Correction Points ▾		Use Master Curve ▾	
	Calibrator	OD	Conc	OD Range	
Point-1	# ▾			Low	High
Point-2	# ▾				
MB Type Factor: ▾		1-Point Calibration Point ▾		<input type="checkbox"/> with Conc-0	
				Stability	
				Reagent Blank # Day # Hour	
				Calibration # Day # Hour	

† = Lot-specific assigned calibrator values.
 # = User defined.

© 2018 Thermo Fisher Scientific, Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AU Series Systems are the registered trademarks of Beckman Coulter.

End