

Infinity Ammonia Application

Beckman Coulter AU680 Parameters

Reagent ID : 154

Specific Test Parameters												
General	LIH	ISE	Range									
Test Name:	Amm		Type:		Plasma ††		Operation:		Yes			
Sample Volume	15		µL	Dilution	0		µL	OD Limit				
Pre-Dilution Rate	1						Min. OD	0.4	Max. O D	2.50		
Reagents Volume	R1 (R1-1)	150		µL	Dilution	0		µL	Reagent OD limit:			
							First	Low	0.8	High	2.50	
							Last	Low	0.8	High	2.50	
	R2 (R2-1)	0		µL	Dilution	0		µL	Dynamic Range Low	10*	High	600*
Common Reagent	Type			Name			Correlation Factor A	1	B	0		
Wavelength:	Pri.	340		Sec.	660		Factor for Maker A	1	B	0		
Method:	FIXED											
Reaction slope:	-						Onboard Stability	14	Days	0	Hour	
Measuring Point 1:	First	2		Last	7							
Measuring Point 2:	First			Last			LIH Influence Check	#				
Linearity:			%				Lipemia	§				
Lag Time Check:							Icterus	§				
							Hemolysis	§				

Specific Test Parameters										
General	LIH	ISE	Range							
Test Name:	AMM		Type:		Plasma ††					
Value/Flag:	VALUE	Level L:	#		Level H:	#				
Specific Ranges:	From		To		Low	High				
	Sex	Year	Month	Year	Month					
1.	#	#	#	#	#	#	#	#	#	#
2.	#	#	#	#	#	#	#	#	#	#
3.	#	#	#	#	#	#	#	#	#	#
4.	#	#	#	#	#	#	#	#	#	#
5.	#	#	#	#	#	#	#	#	#	#
6.	#	#	#	#	#	#	#	#	#	#
7.	No demographics					#	#			
8.	Not within expected values					#	#			
Unit	µmol/L*		Decimal Places	#						

Panic Value

Low	High
#	#

Continued on next page

Infinity Ammonia Application

Beckman Coulter AU680 Parameters

Reagent ID : 154

Calibration Specific									
General									
Test Name:		AMM	Type:		Plasma ††		<input type="checkbox"/> Use Cal.		
Calibration Type:		AB	Formula:		Y=AX+B		Counts:	#	
<Calibrator Parameters>							Slope Check	None	
			Factor Range						
	Calibrator	OD	Conc	Low	High				
Point 1:	#		59 † *	-9999*	9999*				
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				Use Master Curve		<input type="checkbox"/> Lot Calibration	
	Calibrator	OD	Conc	OD Range		Stability			
				Low	High				
Point 1:						Reagent Blanks	7	Day	0
Point 2:						Calibration	7	Day	0
MB Type Factor:		1-Point Calibration Point				<input type="checkbox"/> With CONC-0			

- # User-defined values
 - * Units are in µmol/L. To work in µg/dL multiply by 1.7
 - † Calibrator included in kit
 - § Refer to product IFU for interference information
- ! Do Not Use Ammonium Heparin Plasma Samples**
 †† Sample is EDTA or Lithium Heparin Plasma only

AU680 Contamination Avoidance Parameters							
No.	PRECEDING TEST NAME	FOLLOWING TEST NAME	REAGENT PROBE CLEANER	WASH COUNT	EFFECTIVE OF WATER CLEANING	SAME USE	
						MIXER	CUVETTE
1.	AMM 154	LDH 026	Water	1	Yes	Yes	Yes
2.	GLDH G82100	AMM 154	Water	0	Yes	No	Yes

Note: Other reagents on the carousel which contain/liberate Ammonia may also contaminate Ammonia. Avoid use of the ammonia containing reagents (e.g. Paracetamol OSR6x202) together with OSR61154 to mitigate against atmospheric ammonia transfer. Contact your local Beckman Coulter representative for further information.



WMDE
 Bergerweg 18
 6085 AT Horn
 The Netherlands