

# Infinity Ammonia Application

## Beckman Coulter AU480 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	
	R2 (R2-1)	0		μL	Dilution	0		μL	High
									2.50
									2.50
Wavelength:	Pri.	340		Sec.	660		Dynamic Range Low	10*	
Method:	FIXED			Correlation Factor A	1		High	600*	
Reaction slope:	-			Factor for Maker A	1		B	0	
Measuring Point 1:	First	2		Last	7		Onboard Stability	14 Days	
Measuring Point 2:	First			Last				0 Hour	
Linearity:			%	LIH Influence Check	#		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:	Sex	Year	From Month	To Year	Month	Low	High
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 AU480 Parameters

### Reagent ID 154

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

AU480 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 R1	LDH 026 R2	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.

Note: When selecting a preceding R1 and a following R1, R2 to R2 is also automatically selected. There is no facility for selecting a preceding R2 and a following R2 on the AU480 instrument.



WMDE  
Bergerweg 18  
6085 AT Horn  
The Netherlands