

# Infinity Ammonia Application

## Beckman Coulter AU5800 Parameters

Reagent ID : 154

Parameters		Specific Test Parameters								
General	LIH	ISE	Calculated Tests		Range					
Test Name:	AMM	<	>	Type:	Plasma	Operation:	Yes			
Sample Volume	Volume	<input type="text" value="15"/>	μL	Dilution	<input type="text" value="0"/>	∇μL	OD Limit			
Pre-Dilution Rate		<input type="text" value="1"/>	∇			Min OD	<input type="text" value="0.4"/>	Max OD	<input type="text" value="2.50"/>	
Reagent Volume	R1(R1-1)	<input type="text" value="150"/>	μL	Dilution	<input type="text" value="0"/>	μL	Reagent OD Limit			
	R1-2	<input type="text"/>	μL	Dilution	<input type="text"/>	μL	First Low	<input type="text" value="0.8"/>	High	<input type="text" value="2.50"/>
	R2(R2-1)	<input type="text" value="0"/>	μL	Dilution	<input type="text" value="0"/>	μL	Last Low	<input type="text" value="0.8"/>	High	<input type="text" value="2.50"/>
Common Reagent	Type	<input type="text" value="None"/>		Name	<input type="text" value="None"/>					
Wavelength:	Pri.	<input type="text" value="340"/>	∇nm	Sec.	<input type="text" value="660"/>	∇nm				
Method:		<input type="text" value="FIXED"/>	∇							
Reaction slope:		<input type="text" value="-"/>	∇	Dynamic Range		Low	<input type="text" value="10*"/>	High	<input type="text" value="600*"/>	
Measuring Point 1	First	<input type="text" value="2"/>		Last	<input type="text" value="7"/>	Correlation Factor	A	<input type="text" value="1"/>	B	<input type="text" value="0"/>
Measuring Point 2	First	<input type="text"/>		Last	<input type="text"/>	Factor for Maker	A	<input type="text" value="1"/>	B	<input type="text" value="0"/>
Linearity:		<input type="text"/>	%							
Lag Time Check		<input type="text"/>	∇	On-board Stability Period	<input type="text" value="14"/>	Day	<input type="text" value="#"/>	Hour		
				LIH Influence Check					<input type="text" value="#"/>	
				Lipemia					<input type="text" value="§"/>	
				Icterus					<input type="text" value="§"/>	
				Hemolysis					<input type="text" value="§"/>	

Parameters		Specific Test Parameters						
General	LIH	ISE	Calculated Tests		Range			
Test Name:	AMM	<	>	Type:	Plasma			
Value/Flag:	<input type="text" value="#"/>							
Level				Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>	
Specific Ranges:								
	From		To		Low	High		
<input type="checkbox"/> 1.	Sex	Year	Month	Year	Month	Low	High	
<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							
	8. Not within expected values							
Panic Value	Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>	Unit	<input type="text" value="μmol/L*"/>	Decimal Places	<input type="text" value="#"/>

# Infinity Ammonia Application

## Beckman Coulter AU5800 Parameters

Reagent ID : 154

Parameters		Calibration Parameters			
Calibrators	Calibration Specific	STAT Table Calibration			
Test Name: <input type="text" value="AMM"/> ▾	<input type="text" value="&lt;"/> ▾	Type <input type="text" value="Plasma"/> ▾		Cuvette <input type="text" value=""/> ▾	
<input type="checkbox"/> Use Serum Cal					
Calibration Type: <input type="text" value="AB"/> ▾		Formula: <input type="text" value="Y=AX+B"/> ▾		Counts: <input type="text" value="#"/> ▾	
<b>&lt; Calibrator Parameters &gt;</b>					
Calibrator	OD	Conc	Factor Range Low High	Slope <input type="text" value="None"/> ▾	
Point-1	<input type="text" value="#"/> ▾	<input type="text" value="59†*"/>	<input type="text" value="-9999*"/> <input type="text" value="9999*"/>	Allowable Range Check <input type="checkbox"/> Reagent Blank <input type="text" value=""/> <input type="checkbox"/> Calibration <input type="text" value=""/>  Advanced Calibration Operation <input type="text" value="#"/> ▾ Interval (RB/ACAL) <input type="text" value=""/> ▾  <input type="checkbox"/> Lot Calibration	
Point-2					
Point-3					
Point-4					
Point-5					
Point-6					
Point-7					
Point-8					
Point-9					
Point-10					
<b>&lt;Point Cal. For Master Curve &gt;</b>		No. of Correction Points <input type="text" value=""/> ▾	Use Master Curve <input type="text" value=""/> ▾		
Calibrator	OD	Conc	OD Range Low High	Stability	
Point-1	<input type="text" value=""/> ▾	<input type="text" value=""/>	<input type="text" value=""/> <input type="text" value=""/>	Reagent Blank <input type="text" value="7"/> Day <input type="text" value="#"/> Hour	
Point-2	<input type="text" value=""/> ▾	<input type="text" value=""/>	<input type="text" value=""/> <input type="text" value=""/>	Calibration <input type="text" value="7"/> Day <input type="text" value="#"/> Hour	
MB Type Factor <input type="text" value=""/>	1-Point Calibration Point <input type="text" value=""/> ▾		<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
  - ‡ Sample is EDTA or Lithium Heparin Plasma only
  - § Refer to product IFU for interference information
- ! Do Not Use Ammonium Heparin Plasma Samples**

AU5800 Contamination Avoidance Parameters							
No.	PRECEDING TEST NAME	FOLLOWING TEST NAME	REAGENT PROBE CLEANER	WASH COUNT	CANCEL	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