

Infinity Ammonia Application

Beckman Coulter AU400, AU400e, AU640, AU640e, AU2700, and AU5400 Parameters

Reagent ID 154

Specific Test Parameters										
General	LIH	ISE	Range							
Test Name:	<input type="text" value="AMM"/>		Type:	<input type="text" value="Plasma††"/>		Operation:	<input type="text" value="Yes"/>			
Sample 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. O D	<input type="text" value="2.50"/>	
Reagents Volume	R1 Volume	<input type="text" value="150"/>	μL	Dilution	<input type="text" value="0"/>	μL	Reagent OD limit:			
						First Low	<input type="text" value="0.8"/>	High	<input type="text" value="2.50"/>	
						Last Low	<input type="text" value="0.8"/>	High	<input type="text" value="2.50"/>	
	R2 Volume	<input type="text" value="0"/>	μL	Dilution	<input type="text" value="0"/>	μL	Dynamic Range Low	<input type="text" value="10*"/>	High	<input type="text" value="600*"/>
							Correlation Factor A	<input type="text" value="1"/>	B	<input type="text" value="0"/>
Wavelength:	Pri.	<input type="text" value="340"/>		Sec.	<input type="text" value="660"/>					
Method:	<input type="text" value="FIXED"/>									
Reaction slope:	<input type="text" value="-"/>									
Measuring Point 1:	First	<input type="text" value="2"/>		Last	<input type="text" value="7"/>	Onboard Stability <input type="text" value="14"/> Days				
Measuring Point 2:	First	<input type="text"/>		Last	<input type="text"/>					
Linearity:	<input type="text"/> %									
No Lag Time:	<input type="text"/>									

Specific Test Parameters									
General	LIH	ISE	Range						
Test Name:	<input type="text" value="AMM"/>		Type:	<input type="text" value="Plasma††"/>					
Value/Flag:	<input type="text" value="VALUE"/>	Level L:	<input type="text" value="#"/>	Level H:	<input type="text" value="#"/>				
Specific Ranges:	Sex	From Year	Month	To Year	Month	Low	High	Panic Value Low High <input type="text" value="#"/> <input type="text" value="#"/>	
1.	#	#	#	#	#	#	#		
2.	#	#	#	#	#	#	#		
3.	#	#	#	#	#	#	#		
4.	#	#	#	#	#	#	#		
5.	#	#	#	#	#	#	#		
6.	#	#	#	#	#	#	#		
7.	None selected					#	#		
8.	Out of Range					#	#		
Unit	<input type="text" value="mmol/L*"/>		Decimal Places	<input type="text" value="#"/>					

Continued on next page

Infinity Ammonia Application
Beckman Coulter AU400, AU400e, AU640, AU640e, AU2700, and
AU5400 Parameters
 Reagent ID 154

Calibration Specific								
General	ISE							
Test Name:	AMM	Type:	Plasmatt	<input type="checkbox"/>				
Calibration Type:	AB	Formula:	Y=AX+B	Counts:	#	Process	CONC	
<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:								
1-Point Calibration Point	<input type="checkbox"/>	<input type="checkbox"/>	With CONC-0	Slope Check	None ▾	Advanced Calibration	## ▾	
MB Type Factor	<input type="text"/>					Calibration Stability	7	Day

User-defined values

* Units are in µmol/L. To work in µg/dL multiply by 1.7

† Calibrator included in kit

! Do Not Use Ammonium Heparin Plasma Samples

†† Sample is EDTA or Lithium Heparin Plasma only

AU640/2700/5400 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

AU400 Contamination Avoidance Parameters							
No.	PRECEDING TEST NAME	FOLLOWING TEST NAME	REAGENT PROBE CLEANER	WASH COUNT	CANCEL	SAME USE	
						MIXER	CUVETTE
1.	AMM 154 R1	LDH 026 R2	Water	1	Yes	Yes	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