



Name: Ammonia, DxC 700 AU Ammonia Application

System Reagent: OSR61154
Test name: AMM1G

Reagent ID: 154

General	LIH	ISE	Calculated Test	Range
Test Name: AMM1G				
Type: Plasma ±				
Operation				
Sample Volume	15 µL	Dilution	0 µL	OD Limit
Pre-Dilution Rate	1			Min. OD 0.4 Max OD 2.50
Reagent Volume R1 (R1-1)	150 µL	Dilution	0 µL	Reagent OD Limit 1 st Low 0.8 High 2.50
R1-2		Dilution		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 nm	Sec	660 nm	Manufacturer Factor A 1 B 0
Method	FIXED			
Reaction Slope	-			Onboard Stability Period 14 Day 0 Hour
Measuring Point-1 1st	2	Last	7	LIH Influence Check #
Measuring Point-2 1st		Last		Lipemia £
Linearity Limit				Icterus £
Lag Time Check				Hemolysis £

General	LIH	ISE	Calculated Test	Range
Test Name: AMM1G				
Type: Plasma ±				
Value Flag	#	Level	Low # High #	
Specific Ranges				
	From	To	Other Type	Low High
Sex	Year Month	Year Month	None	# #
□ 1:	# #	# #	None	# #
□ 2:	# #	# #	None	# #
□ 3:	# #	# #	None	# #
□ 4:	# #	# #	None	# #
□ 5:	# #	# #	None	# #
□ 6:	# #	# #	None	# #
7:	Standard demographics			# #
8:	Not within expected values			# #
Panic Value	Low #	High #	Unit µmol/L*	Select Decimal Places #

Calibrators	General	ISE
Test Name: AMM1G		
Type: Plasma±		
<input type="checkbox"/> Use Serum Cal.		
Calibration Type:	AB	Formula: Y=AX+B
		Counts: #
<Calibrator Parameters>		Slope Check None
	Calibrator	OD
		Conc
		Range
		Low High
Point-1	#	
Point-2		59 * ±±
Point-3		
Point-4		
Point-5		
Point-6		
Point-7		
MB Type Factor		1-Point Calibration Point
		<input type="checkbox"/> with Conc-0
		Stability
		Reagent Blank 7 Day 0 Hour
		Calibration 7 Day 0 Hour
		Interval (RB)
		Interval (ACAL)
		Allowable Range Check
		<input type="checkbox"/> Reagent Blank #
		<input type="checkbox"/> Calibration #
		Advanced Calibration
		Operation #

User Defined

* Values set for working in µmol/L. To work in µg/dL multiply by 1.7

C - Closed, Value/Text must be defined in the setting sheet

P - Pre-populate, editable, Value/Text must be defined in the setting sheet

±- Sample is only EDTA plasma or Heparin Plasma. Do not use Ammonia Heparin Plasma Samples.

£- Refer to IFU for interference information.

±±- Calibrator included in kit

DxC 700 AU Contamination Prevention Parameters							
No.	PRECEDING TEST NAME	FOLLOWING TEST NAME	REAGENT PROBE CLEANER	WASH COUNT	CANCEL	PREVENT 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.