

Infinity™ Ammonia Application

Beckman Coulter AU480 Parameters

Reagent ID 154

Catalog No. **REF** 106518This Application is Intended for the in vitro quantitative determination of Ammonia (NH₃) in Human Plasma**IVD**

For In Vitro Diagnostic Use Only

Rx Only

The information provided in this application sheet is intended as a supplement to the product Instructions for Use (IFU). Refer to the package IFU on intended use, reagent storage, reagent preparation, specimen collection, specimen preparation, specimen storage, quality control, and additional performance data.



**Fisher Diagnostics, a division of Fisher Scientific Company, LLC,
a part of Thermo Fisher Scientific Inc,
8365 Valley Pike, Middletown, VA 22645-1905 USA**



B-R-A-H-M-S GmbH, Neuendorfstrasse 25, 16761 Hennigsdorf, Germany

**Produced by Fisher Diagnostics for:
Beckman Coulter Inc.
250 S. Kraemer Blvd.
Brea, CA 92821 USA**

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		
Reagents Volume	R1 (R1-1)	150		µL	Dilution	0		µL	Max. O D	2.50	
	R2 (R2-1)	0		µL	Dilution	0		µL	Reagent OD limit:		
Wavelength:	Pri.	340			Sec.	660			First Low	0.8	
	Method:	FIXED							Last Low	0.8	
Reaction slope:	-								High	2.50	
Measuring Point 1:	First	2			Last	7			High	2.50	
Measuring Point 2:	First				Last				Dynamic Range Low	10*	
Linearity:			%						Correlation Factor A	1	
Lag Time Check:									Factor for Maker A	1	
									High	600*	
									Onboard Stability	14 Days	
										0 Hour	
									LIH Influence Check	#	
									Lipemia	§	
									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	Month	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

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:	<input type="text" value="#"/>		<input type="text" value="59 † *"/>	<input type="text" value="-9999*"/>	<input type="text" value="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		Use Master Curve	<input type="text"/>		<input type="checkbox"/> Lot Calibration		
				Low	High	Stability					
Point 1:						Reagent Blanks	<input type="text" value="7"/>	Day	<input type="text" value="0"/>		
Point 2:						Calibration	<input type="text" value="7"/>	Day	<input type="text" value="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

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 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.