

# Infinity™ Ammonia Application

## Beckman Coulter DxC 700 AU Analyzer

Catalog No. **REF** 106518

This Application is Intended for the in vitro quantitative determination of Ammonia (NH<sub>3</sub>) 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**

W P	W P	Test No	Test No	Ser/Uri/Ot	Ser/Uri/Ot	Yes
-----	-----	---------	---------	------------	------------	-----

### Name: Ammonia, DxC 700 AU Ammonia Application

System Reagent: OSR61154  
 Test name: AMM1G

Reagent ID: 154

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

Calibrators	General	ISE
<b>Test Name:</b> AMM1G		
<b>Type:</b> 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

# User Defined \* Values set for working in µmol/L. To work in µg/dL multiply by 1.7 £- Refer to IFU for interference information.  
 C - Closed, Value/Text must be defined in the setting sheet ±±- Calibrator included in kit  
 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.**

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 together with OSR61154 to mitigate against atmospheric ammonia transfer. Contact your local Beckman Coulter representative for further information.