Infinity™ AST (GOT) Liquid Stable Reagent

PRODUCT SUMMARY

- Stability Linear Range Specimen Type Method Reagent Preparation :
- Until Expirv at 2-8°C Up to 450 U/L (7.52 µkat/L) Serum Kinetic Supplied ready to use.

IVD INTENDED USE

This reagent is intended for the in vitro quantitative determination of AST (Aspartate Aminotransferase EC2.6.1.1) in human serum.

CLINICAL SIGNIFICANCE

AST is widely distributed with high concentrations in the heart, liver, skeletal muscle, kidney and erythrocytes. Damage or disease to any of these tissues such as myocardial infarction, viral hepatitis, liver necrosis, cirrhosis and muscular dystrophy may result in raised serum levels of AST.¹

METHODOLOGY

In 1955, Karmen et al² described the first kinetic assay of AST for diagnostic purposes. This method was evaluated and improved by many investigators primarily Henry et al³ and now forms the basis of many national and international recommended procedures. This AST Reagent is based on the recommendations of the IFCC.4

The series of reactions involved in the assay system is as follows:

1. L-Aspartate + 2-Oxoglutarate \xrightarrow{AST} Oxaloacetate + L-Glutamate

- 3. Sample Pyruvate + NADH _____ L-Lactate + NAD
- AST present in the sample catalyses the transfer of the amino group from 1 L-aspartate to 2-oxoglutarate forming oxaloacetate and L-glutamate.
- Oxaloacetate in the presence of NADH and Malate dehydrogenase (MDH), is 2 reduced to L-malate. In this reaction NADH is oxidized to NAD. The reaction is monitored by measuring the rate of decrease in absorbance at 340nm due to the oxidation of NADH to NAD.
- Addition of Lactate dehydrogenase (LDH) to the reagent is necessary to 3 achieve rapid and complete reduction of endogenous pyruvate so that it does not interfere with the assay.

REAGENT COMPOSITION

Active Ingredients	Concentration
2-Oxoglutarate	13 mmol/L
L-Aspartate	220 mmol/L
MDH (microbial)	> 100 U/L
LDH (microbial)	> 1500 U/L
NADH	> 0.12 mmol/L
Tris Buffer	88 mmol/L
EDTA	5.0 mmol/L

pH 8.10 ± 0.1 at 20°C.

WARNING: Do not ingest. Avoid contact with skin and eyes. If spilt, thoroughly wash affected areas with water. Reagent contains Sodium Azide which may react with copper or lead plumbing. Flush with plenty of water when disposing. For further information consult the Infinity AST(GOT) Liquid Stable Reagent Material Safety Data Sheet.

REAGENT PREPARATION

Reagent is supplied ready to use.

SYMBOLS IN PRODUCT LABELLING

- EC REP Authorized Representative For in vitro diagnostic use IVD Batch code/Lot number LOT REF Catalogue number
- X **Temperature Limitation**
 - Use by/Expiration Date

Manufactured by

CAUTION. CONSULT INSTRUCTIONS FOR USE.

i Consult instructions for use

STABILITY AND STORAGE

When stored at 2-8°C reagent is stable until the expiration date stated on the bottle and kit box label. It is recommended that when the reagent is not in use for prolonged periods of time (eg: overnight) the reagent be capped and stored at 2-8°C. Indications of Reagent Deterioration:

- Turbidity,
- Absorbance < 1.0 at 340 nm (1 cm),
- Failure to recover control values within the assigned range.

SPECIMEN COLLECTION AND HANDLING

Use non-haemolysed serum. Serum

AST samples may be stored for at least 7 days at 4°C.5 Storage:

ADDITIONAL EQUIPMENT REQUIRED BUT NOT PROVIDED

- If required, pipettes for accurately dispensing measured volumes. A clinical chemistry analyzer capable of maintaining constant temperature (37°C)
- and measuring absorbance at 340 nm.
- Analyzer specific consumables, eg: sample cups. Normal and Abnormal control material.

ASSAY PROCEDURE

The following system parameters are recommended. Individual instrument applications are available upon request from the Technical Support Group.

SYSTEM PARAMETERS

Temperat Primary V Seconda Assay Ty	ture Vavelength ry Wavelength pe	37°C 340 nm 405 nm Bate/Kii	(334, 365nm) netic
Direction		Decreas	se
Sample :	Reagent Ratio	1:10	
eg:	Sample Vol	30 µL	
	Reagent Vol	300 µL	
Delay/Lag	g	30 seco	nds
Read Tirr	ne	1 to 3 m	ninutes
Reagent	Blank	Low	1.0 AU
(1cm ligh	tpath, 340nm)	High	2.5 AU
Linearity		450 U/L	. (7.52 µkat/L)
(refer to L	inearity section)		
Analytica	I Sensitivity	0.573 ∆	mA/min per U/L
(1cm ligh	tpath, 340nm)	(34.31 /	ΔmA/min per μkat/L)

CALCULATIONS

Results are calculated, usually automatically by the instrument, as follows:

Activity in U/L = \triangle Abs/min x Factor

Factor =
$$\frac{\text{TV x 1000}}{6.3 \text{ x SV x P}}$$

Where: TV

- Total reaction volume in mL = Sample volume in mL
- SV
- millimolar absorption coefficent of NADH at 340nm (See note 4). 6.3 Р
- Cuvette pathlength in cm.



Example:

 $\Delta Abs/min = 0.08$

= 1746 Factor = 0.08 x 1746 = 140 U/L AST

NOTES

- 1. The reagent and sample volumes may be altered proportionally to accommodate different spectrophotometer requirements.
- 2. If the change in absorbance is greater than 0.26/min repeat the assay with less sample or dilute with saline. Remember to adjust the factor for the smaller sample volume or to multiply the final result by the dilution factor.
- 3. Valid results depend on an accurately calibrated instrument, timing, and temperature control.
- The millimolar absorption coefficent for NADH at 334nm = 6.18 and at 365nm 4. = 3.40
- 5. Unit Conversion: U/L x 16.67 x 10⁻³ = µkat/L

CALIBRATION

Not required. The rate of reaction is converted to U/L of activity by a calculation factor. Refer to the calculation section of this package insert.

QUALITY CONTROL

To ensure adequate quality control, normal and abnormal control with assayed values should be run as unknown samples:-

- At least once per day or as established by the laboratory.
- When a new bottle of reagent is used.
- After preventative maintenance is performed or a critical component is . replaced

Control results falling outside the upper or lower limits of the established ranges indicate the assay may be out of control.

- The following corrective actions are recommended in such situations:-
- Repeat the same controls. If repeated control results are outside the limits, prepare fresh control serum
- and repeat the test. If results on fresh control material still remain outside the limits, then repeat the test with fresh reagent.
- If results are still out of control, contact Technical Services or the local distributor.

LIMITATIONS

Studies to determine the level of interference from bilirubin (free & conjugated), 1. haemoglobin and lipaemia were carried out using commercially available interference check products. The following results were obtained:

Haemoglobin: No interference from haemoglobin up to a level of 150 mg/dL.

Free bilirubin: No interference from free bilirubin up to a level of 260 umol/L (15 mg/dL).

Conjugated bilirubin: No interference from conjugated bilirubin up to a level of 116 µmol/L (6.8 mg/dL).

Lipaemia: No interference from lipaemia, measured as an absorbance at 630nm, up to 1.68 AU.

- Haemolyzed serum specimens should not be used. AST activity levels in 2. erythrocytes are some 15 times higher, than those in sera.6
- Young DS7 has published a comprehensive list of drugs and substances З. which may interfere with this assay.

EXPECTED VALUES⁵

At 37°C 5 - 34 U/L (0.084 - 0.568 µkat/L)

Levels approximately twice the adult levels are seen in neonates and infants. These levels decline to normal adult levels after 6 months.

The quoted values should serve as a guide only. It is recommended that each laboratory verify this range or derives a reference interval for the population that it serves.8

PERFORMANCE DATA

The following data was obtained using the Infinity AST(GOT) Liquid Stable Reagent on a well maintained automated clinical chemistry analyzer. Users should establish product performance on their specific analyzer used.

IMPRECISION

Imprecision was evaluated using two levels of commercial control and following the NCCLS EP5-T procedure⁹.

		LEVEL I	LEVEL II
Number of	data points	80	80
Mean (U/L / µkat/L)		45 / 0.752	194 / 3.24
Within Run: SD (U/L / µkat/L)		0.7 / 0.012	1.4 / 0.023
	CV (%)	1.6	0.7
Total:	SD(U/L / µkat/L)	1.2 / 0.020	2.3 / 0.038
	CV (%)	2.7	1.2

METHOD COMPARISON

Comparison studies were carried out using a similar commercially available reagent as a reference. Serum samples were assayed in parallel and the results compared by least squares regression. The following statistics were obtained.

Number of sample pairs	84
Range of sample results	8 - 276 U/L(0.134 - 4.61 µkat/L)
lean of reference method results	37 U/L (0.618 µkat/L)
Aean of Infinity AST (GOT) results	38 U/L (0.635 µkat/L)
Slope	0.98
ntercept	2.1 U/L (0.035 µkat/L)
Correlation coefficient	0.997

LINEARITY

When run as recommended, the assay is linear up to 450 U/L (7.52 $\mu kat/L).$ Linearity on automated instruments will be dependent upon the ratio of sample volume to reagent volume used and the timing of measurements. The specific instrument application should be consulted.

ANALYTICAL SENSITIVITY

When run as recommended the sensitivity of this assay is 0.573 Δ mA /min per U/L (34.31 AmA /min per µkat/L).

REFERENCES

- Zilva JF, Pannall PR. "Plasma Enzymes in Diagnosis" in Clinical Chemistry in 1. Diagnosis and Treatment. Lloyd-Luke London. 1979; Chap 15:338-9.
- 2 Karmen A. J Clin Investigation 1955; 43:131.
- Henry RJ, et al. Am J Clin Path 1960; 34:381.
- 4. IFCC Method for L-Aspartate aminotransferase. J. Clin. Chem. Clin. Biochem. 1986: 24:497-510.
- Murray RL. "Aspartate aminotransferase" in Clinical Chemistry. Theory, 5 analysis and correlation. Kaplan LA, Pesce AJ (Ed), CV Mosby Company 1984; 1105-8.
- Burtis CA, Ashwood ER, "Tietz textbook of Clinical Chemistry" Second Edition, 6. 1994; 795.
- 7 Young DS, Effects of Drugs on Clinical Laboratory Tests. Third Edition. 1990: 3.45-52
- 8. Wachtel M et al, Creation and Verification of Reference Intervals. Laboratory Medicine 1995; 26:593-7.
- National Committee for Clinical Laboratory Standards. User evaluation of Precision Performance of Clinical Chemistry devices. NCCLS; 1984, NCCLS Publication EP5-T

Fisher Diagnostics a division of Fisher Scientific Company, LLC a part of Thermo Fisher Scientific Inc. Middletown, VA 22645-1905 USA Phone: 800-528-0494 540-869-3200 Fax: 540-869-8132

MDCI Ltd. ECREP Arundel House 1 Liverpool Gardens Worthing, West Sussex BN11 1SL UK © 2010 Thermo Fisher Scientific Inc. All rights reserved.

REF

Reorder Information **Configuration**

TR70121

2 x 125 mL

Catalogue No.