



**IMMUNOASSAYS AND SERVICES**

**BIOGENIC AMINES & NEUROSCIENCE | ENDOCRINOLOGY | FOOD SAFETY**

LABOR DIAGNOSTIKA NORD GmbH & Co.KG | Am Eichenhain 1 | 48531 Nordhorn | Germany | Tel. +49 5921 8197-0 | Fax +49 5921 8197-222 | info@ldn.de | www.ldn.de

**Instructions for use**  
**Tryptophan ELISA**  
**Assay Kit**

Distributed in the US and Canada by:

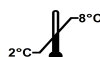
**EAGLE BIOSCIENCES, INC.**  
20A NW BLVD, SUITE 112 NASHUA, NH 03063  
P: 617-419-2019 F: 617-419-1110  
WWW.EAGLEBIO.COM — INFO@EAGLEBIO.COM



**EAGLE**  
BIOSCIENCES

**REF**

**BA E-2700**



**RUO**

**CE**

## 1. Introduction

### 1.1 Intended use and principle of the test

Enzyme Immunoassay for the quantitative determination of Tryptophan in urine, serum and plasma samples.

After extraction and derivatization Tryptophan is quantitatively determined by ELISA.

The competitive ELISA uses the microtiter plate format. The antigen is bound to the solid phase of the microtiter plate. The derivatized standards, controls and samples and the solid phase bound analyte compete for a fixed number of antibody binding sites. When the system is in equilibrium, free antigen and free antigen-antibody complexes are removed by washing. The antibody bound to the solid phase is detected by an anti-rabbit IgG-peroxidase conjugate using TMB as a substrate. The reaction is monitored at 450 nm.

Quantification of unknown samples is achieved by comparing their absorbance with a standard curve prepared with known standards.

### 1.2 Clinical application

L-Tryptophan is one of the essential amino acids for the human metabolism and must be part of its diet.

In humans it serves as precursor for the synthesis of the neurotransmitters serotonin and tryptamine as well as for the synthesis of nicotinic acid and the epiphyseal hormone melatonin. Tryptophan is catabolized to kynurenine through the enzyme IDO (indoleamine-2,3-dioxygenase). Increased IDO activity is an expression of neuro-endocrine-immunological dysregulation, which is often associated with depressive symptoms such as bipolar disorder (manic depression). In addition Tryptophan and its metabolites regulate neurobehavioral effects such as appetite, sleeping-waking-rhythm and pain perception.

This kit is for research use only and should not be used as a diagnostic tool.

## 2. Procedural cautions, guidelines, warnings and limitations

### 2.1 Procedural cautions, guidelines and warnings

- (1) This kit is intended for professional use only. Users should have a thorough understanding of this protocol for the successful use of this kit. Only the test instruction provided with the kit is valid and has to be used to run the assay. Reliable performance will only be attained by strict and careful adherence to the instructions provided.
- (2) This assay was validated for certain types of samples as indicated in *Intended Use* (please refer to Chapter 1). Any off-label use of this kit is in the responsibility of the user and the manufacturer cannot be held liable.
- (3) The principles of Good Laboratory Practice (GLP) have to be followed.
- (4) In order to reduce exposure to potentially harmful substances, wear lab coats, disposable protective gloves and protective glasses where necessary.
- (5) All kit reagents and specimens should be brought to room temperature and mixed gently but thoroughly before use. Avoid repeated freezing and thawing of reagents and specimens.
- (6) For dilution or reconstitution purposes, use deionized, distilled or ultra-pure water.
- (7) The microplate contains snap-off strips. Unused wells must be stored at 2 °C to 8 °C in the sealed foil pouch with desiccant and used in the frame provided.
- (8) Duplicate determination of sample is highly recommended to be able to identify potential pipetting errors.
- (9) Once the test has been started, all steps should be completed without interruption. Make sure that the required reagents, materials and devices are prepared ready at the appropriate time.
- (10) Incubation times do influence the results. All wells should be handled in the same order and time intervals.
- (11) To avoid cross-contamination of reagents, use new disposable pipette tips for dispensing each reagent, sample, standard and control.
- (12) A standard curve must be established for each run.
- (13) The controls should be included in each run and fall within established confidence limits. The confidence limits are listed in the QC-Report.
- (14) Do not mix kit components with different lot numbers within a test and do not use reagents beyond expiry date as shown on the kit labels.
- (15) Avoid contact with Stop Solution containing 0.25 M H<sub>2</sub>SO<sub>4</sub>. It may cause skin irritation and burns. In case of contact with eyes or skin, rinse off immediately with water.

- (16) TMB substrate has an irritant effect on skin and mucosa. In case of possible contact, wash eyes with an abundant volume of water and skin with soap and abundant water. Wash contaminated objects before reusing them.
- (17) For information on hazardous substances included in the kit please refer to Safety Data Sheet (SDS). The Safety Data Sheet for this product is made available directly on the website of the manufacturer or upon request.
- (18) The expected reference values reported in this test instruction are only indicative. It is recommended that each laboratory establishes its own reference intervals.
- (19) This kit is for research use only and should not be used as a diagnostic tool.
- (20) Kit reagents must be regarded as hazardous waste and disposed of according to national regulations.

## 2.2 Limitations

Any inappropriate handling of samples or modification of this test might influence the results.

### 2.2.1 Interfering substances

#### Serum/Plasma

Samples containing precipitates or fibrin strands or which are haemolytic or lipemic might cause inaccurate results.

#### 24-hour urine

Please note the sample preparation! If the percentage of the final concentration of acid is too high, this will lead to incorrect results for the urine samples.

### 2.2.2 Drug interferences

There are no known substances (drugs) which ingestion interferes with the measurement of tryptophan level in the sample.

### 2.2.3 High-Dose-Hook effect

No hook effect was observed in this test.

## 3. Storage and stability

Store the unopened reagents at 2 - 8 °C until expiration date. Do not use components beyond the expiry date indicated on the kit labels. Once opened the reagents are stable for 1 month when stored at 2 - 8 °C. Once the resealable pouch has been opened, care should be taken to close it tightly with desiccant again.


## 4. Materials

### 4.1 Contents of the kit

<b>BA D-0090</b>	<b>FOILS</b>	<b>Adhesive Foil</b> - Ready to use
Contents:	Adhesive Foils in a resealable pouch	
Volume:	1 x 4 foils	
<b>BA D-0024</b>	<b>REAC-PLATE</b>	<b>Reaction Plate</b> - Ready to use
Contents:	1 x 96 well plate, empty in a resealable pouch	
<b>BA E-0030</b>	<b>WASH-CONC 50x</b>	<b>Wash Buffer Concentrate</b> - Concentrated 50x
Contents:	Buffer with a non-ionic detergent and physiological pH	
Volume:	1 x 20 ml/vial, light purple cap	
<b>BA E-0040</b>	<b>CONJUGATE</b>	<b>Enzyme Conjugate</b> - Ready to use
Contents:	Goat anti-rabbit immunoglobulins conjugated with peroxidase	
Volume:	1 x 12 ml/vial, red cap	
<b>BA E-0055</b>	<b>SUBSTRATE</b>	<b>Substrate</b> - Ready to use
Contents:	Chromogenic substrate containing tetramethylbenzidine, substrate buffer and hydrogen peroxide	
Volume:	1 x 12 ml/black vial, black cap	

**BA E-0080** **STOP-SOLN** **Stop Solution** - Ready to use

Contents: 0.25 M sulfuric acid  
 Volume: 1 x 12 ml/vial, light grey cap

Hazards identification: 

H290 May be corrosive to metals.  
 H314 Causes severe skin burns and eye damage.

**BA E-2731** **TRYP** **Tryptophan Microtiter Strips** - Ready to use

Contents: 1 x 96 well (12x8) antigen precoated microwell plate in a resealable pouch with desiccant

**BA E-2710** **AS TRYP** **Tryptophan Antiserum** - Ready to use

Contents: Rabbit anti-tryptophan antibody, blue coloured  
 Volume: 1 x 6 ml/vial, blue cap

**Standards and Controls** - Ready to use

Cat. no.	Component	Colour/Cap	Concentration µg/ml	Concentration µmol/l	Volume/ Vial
<b>BA E-2701</b>	<b>STANDARD A</b>	white	0	0	4 ml
<b>BA E-2702</b>	<b>STANDARD B</b>	light yellow	2.5	12.2	4 ml
<b>BA E-2703</b>	<b>STANDARD C</b>	orange	7.5	36.7	4 ml
<b>BA E-2704</b>	<b>STANDARD D</b>	dark blue	25	122	4 ml
<b>BA E-2705</b>	<b>STANDARD E</b>	light grey	75	367	4 ml
<b>BA E-2706</b>	<b>STANDARD F</b>	black	250	1 224	4 ml
<b>BA E-2751</b>	<b>CONTROL 1</b>	light green	Refer to QC-Report for expected value and acceptable range!		4 ml
<b>BA E-2752</b>	<b>CONTROL 2</b>	dark red			4 ml

Conversion: Tryptophan (µg/ml) x 4.89 = Tryptophan (µmol/l)

Contents: Acidic buffer with non-mercury stabilizer, spiked with defined quantity of tryptophan

**BA E-2413** **ASSAY-BUFF** **Assay Buffer** - Ready to use

Contents: Buffer with alkaline pH  
 Volume: 1 x 20 ml/vial, yellow cap

**BA E-2428** **EQUA-REAG** **Equalizing Reagent** - Lyophilized

Contents: Lyophilized protein  
 Volume: 1 vial, brown cap

**BA E-2446** **D-REAGENT** **D-Reagent** - Ready to use

Contents: Crosslinking agent in dimethylsulfoxide  
 Volume: 1 x 4 ml/vial, white cap

**BA E-2458** **Q-BUFFER** **Q-Buffer** - Ready to use

Contents: TRIS buffer  
 Volume: 1 x 20 ml/vial, white cap

**BA E-2788** **PBS** **PBS** - Ready to use

Contents: Phosphate Buffered Saline  
 Volume: 1 x 20 ml/vial, orange cap

**BA E-2721** **PREC-REAG** **Precipitating Reagent** - Ready to use

Contents: Acidic reagent for precipitation of plasma/serum proteins, red coloured  
 Volume: 1 x 4 ml/vial, white cap

#### 4.2 Additional materials and equipment required but not provided in the kit

- Calibrated precision pipettes to dispense volumes between 10 – 300 µl; 12.5 ml
- Polystyrene or polypropylene tubes and suitable rack
- Microtiter plate washing device (manual, semi-automated or automated)
- ELISA reader capable of reading absorbance at 450 nm and if possible 620 - 650 nm
- Microtiter plate shaker (shaking amplitude 3 mm; approx. 600 rpm)
- Absorbent material (paper towel)
- Water (deionized, distilled or ultra-pure)
- Vortex mixer

#### 5. Sample collection and storage

##### **Plasma**

Whole blood should be collected by venipuncture into centrifuge tubes containing EDTA as anti-coagulant (Monovette™ or Vacuette™ for plasma) and centrifuged according to manufacturer's instructions at room temperature immediately after collection.

Fasting specimens or pre-feed specimens for children (2 - 3 hours after last meal) are advised.

Haemolytic and especially lipemic samples should not be used for the assay.

Storage: up to 48 hours at 2 - 8 °C, for longer period (up to 6 month) at -20 °C.

Repeated freezing and thawing should be avoided.

##### **Serum**

Collect blood by venipuncture (Monovette™ or Vacuette™ for serum), allow to clot, and separate serum by centrifugation according to manufacturer's instructions at room temperature. Do not centrifuge before complete clotting has occurred. Patients receiving anticoagulant therapy may require increased clotting time.

Fasting specimens or pre-feed specimens for children (2 - 3 hours after last meal) are advised.

Haemolytic and especially lipemic samples should not be used for the assay.

Storage: up to 48 hours at 2 - 8 °C, for longer period (up to 6 month) at -20 °C.

Repeated freezing and thawing should be avoided.

##### **Urine**

Spontaneous urine or 24-hour urine, collected in a bottle containing 10 - 15 ml of 6 M HCl, can be used.

*If 24-hour urine is used please record the total volume of the collected urine.*

Storage: for longer periods (up to 6 month) at -20 °C.

Repeated freezing and thawing should be avoided. Avoid exposure to direct sunlight.

#### 6. Test procedure

Allow all reagents and samples to reach room temperature and mix thoroughly by gentle inversion before use. Duplicate determinations are recommended. It is recommended to number the strips of the microwell plate before usage to avoid any mix-up.

The binding of the antisera and of the enzyme conjugate and the activity of the enzyme are temperature dependent, and the absorbance values may vary if a thermostat is not used. The higher the temperature, the higher the extinction values will be. Corresponding variations also apply to the incubation times. The optimal temperature during the Enzyme Immunoassay is between 20 – 25 °C.

#### 6.1 Preparation of reagents

##### **Wash Buffer**

Dilute the 20 ml Wash Buffer Concentrate with water (deionized, distilled or ultra-pure) to a final volume of 1000 ml.

Storage: 1 month at 2 – 8 °C

##### **Equalizing Reagent**

Reconstitute the Equalizing Reagent with **12.5 ml** of **Assay Buffer**.

Reconstituted Equalizing Reagent which is not used immediately has to be stored in aliquots for max 1 month at -20 °C and may be thawed only once.


##### **D-Reagent**

The D-Reagent has a freezing point of 18.5 °C. It must be ensured that the D-Reagent has reached room temperature and forms a homogeneous, crystal-free solution.

##### **Tryptophan Microtiter Strips**

In rare cases residues of the blocking and stabilizing reagent can be seen in the wells as small, white dots or lines. These residues do not influence the quality of the product.

## 6.2 Precipitation

1.	Pipette <b>20 µl</b> of the <b>standards, controls</b> and <b>samples</b> into the respective <b>tubes</b> .
2.	Add <b>200 µl PBS</b> to all tubes.
3.	Add <b>25 µl Precipitating Reagent</b> to all tubes.
4.	Mix the <b>tubes</b> thoroughly (vortex) and centrifuge for <b>15 minutes</b> at <b>3,000 x g</b> .
	Take <b>25 µl</b> of the clear supernatant for the <b>derivatization</b> .

## 6.3 Derivatization

1.	Pipette <b>25 µl</b> of the <b>precipitated standards, controls</b> and <b>samples</b> into the appropriate wells of the <b>Reaction Plate</b> .
2.	Pipette <b>50 µl</b> of the <b>Equalizing Reagent</b> into all wells.
3.	Pipette <b>10 µl</b> of the <b>D-Reagent</b> into all wells.
4.	Cover plate with <b>Adhesive Foil</b> and incubate for <b>2 h</b> at <b>RT</b> (20 – 25 °C) on a <b>shaker</b> (approx. 600 rpm).
5.	Pipette <b>100 µl</b> of the <b>Q-Buffer</b> into all wells.
6.	Incubate for <b>10 min</b> at <b>RT</b> (20 – 25 °C) on a <b>shaker</b> (approx. 600 rpm).
7.	<b>Use 25 µl for the ELISA!</b>


## 6.4 Tryptophan ELISA

1.	Pipette <b>25 µl</b> of the <b>prepared standards, controls and samples</b> into the appropriate wells of the <b>Tryptophan Microtiter Strips</b> .
2.	Pipette <b>50 µl</b> of the <b>Tryptophan Antiserum</b> into all wells and mix shortly.
3.	Cover plate with <b>Adhesive Foil</b> and incubate for <b>15 - 20 h</b> (overnight) at <b>2 – 8 °C</b> .
4.	Remove the foil. Discard or aspirate the content of the wells. Wash the plate <b>3 x</b> by adding <b>300 µl</b> of <b>Wash Buffer</b> , <b>discarding</b> the content and <b>blotting dry each time</b> by tapping the inverted plate on absorbent material.
5.	Pipette <b>100 µl</b> of the <b>Enzyme Conjugate</b> into all wells.
6.	Incubate for <b>30 min</b> at <b>RT</b> (20 – 25 °C) on a <b>shaker</b> (approx. 600 rpm).
7.	Discard or aspirate the content of the wells. Wash the plate <b>3 x</b> by adding <b>300 µl</b> of <b>Wash Buffer</b> , <b>discarding</b> the content and <b>blotting dry each time</b> by tapping the inverted plate on absorbent material.
8.	Pipette <b>100 µl</b> of the <b>Substrate</b> into all wells and incubate for <b>20 - 30 min</b> at <b>RT</b> (20 – 25 °C) on a <b>shaker</b> (approx. 600 rpm). <b>Avoid exposure to direct sunlight!</b>
9.	Add <b>100 µl</b> of the <b>Stop Solution</b> to each well and shake the microtiter plate to ensure a homogeneous distribution of the solution.
10.	<b>Read</b> the <b>absorbance</b> of the solution in the wells within 10 minutes, using a microplate reader set to <b>450 nm</b> (if available a reference wavelength between 620 nm and 650 nm is recommended).

## 7. Calculation of results

<b>Measuring range</b>	<b>Tryptophan</b>
	1.2 - 250 µg/ml

The calibration curve is obtained by plotting the absorbance readings (calculate the mean absorbance) of the standards (linear, y-axis) against the corresponding standard concentrations (logarithmic, x-axis).  
Use non-linear regression for curve fitting (e.g. spline, 4-parameter, akima).

 *This assay is a competitive assay. This means: the OD-values are decreasing with increasing concentrations of the analyte. OD-values found below the standard curve correspond to high concentrations of the analyte in the sample and have to be reported as being positive.*

The concentrations of the samples and controls can be read directly from the standard curve.

The total amount of Tryptophan excreted in urine during 24 h is calculated as following:

$$\mu\text{g}/24\text{h} = \mu\text{g}/\text{l} \times \text{l}/24\text{h}$$

### Conversion

$$\text{Tryptophan } (\mu\text{g}/\text{ml}) \times 4.89 = \text{Tryptophan } (\mu\text{mol}/\text{l})$$

### Expected reference values


It is strongly recommended that each laboratory should determine its own reference values.

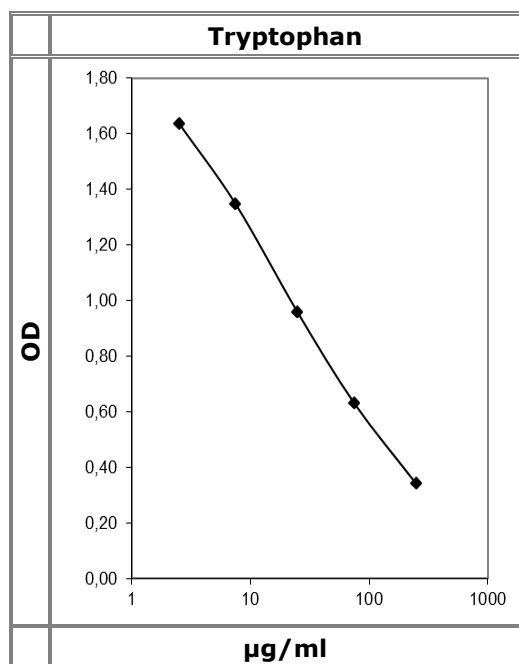
<b>Plasma / Serum</b>	<b>Spontaneous urine</b>
9.3 - 17 µg/ml	1.5 - 40 µg/g creatinine

### 7.1 Quality control

It is recommended to use control samples according to national regulations. Use controls at both normal and pathological levels. The kit controls or other commercial controls should fall within established confidence limits. The confidence limits of the kit controls are indicated on the QC-Report.

### 7.2 Typical standard curve

 *Example, do not use for calculation!*



## 8. Assay characteristics

<b>Analytical Sensitivity (Limit of Detection)</b>	<b>Tryptophan</b>
	1.2 µg/ml

<b>Analytical Specificity (Cross Reactivity)</b>	<b>Substance</b>	<b>Cross Reactivity (%)</b>
	Tryptophan	100
	5-Hydroxy-L-tryptophan	<0.01
	Tryptamine	<0.01
	5-Methoxy-L-tryptophan	<0.01
	5-Hydroxytryptamine	<0.01
	5-Methoxytryptamine	<0.01

<b>Precision</b>					
<b>Intra-Assay</b>			<b>Inter-Assay</b>		
Sample	Range (µg/ml)	CV (%)	Sample	Range (µg/ml)	CV (%)
1 (n = 77)	9.4 ± 1.0	11	1 (n = 16)	9.2 ± 1.4	15
2 (n = 78)	27 ± 2.8	11	2 (n = 16)	45 ± 4	8.4

<b>Linearity</b>		Range (µg/ml)	Serial dilution up to	Range (%)
	Urine		1.3 - 100	1:75







<b>Recovery</b>		Mean (%)	Range (%)
	Urine	106	104 - 110
	Serum	95	86 - 100

## 9. References/Literature

- (1) El-Bakly et al. Hypericum Perforatum Decreased Hippocampus TNF-α and Corticosterone Levels with No Effect on Kynurenine/Tryptophan Ratio in Bilateral Ovariectomized Rats. Korean J Physiol Pharmacol, 18:133-139 (2014)
- (2) Nowak et al. Tryptophan hydroxylase-1 regulates immune tolerance and inflammation. The Journal of Experimental Medicine, 209(11): 2127-2135 (2012)
- (3) Sorensen et al. Indoleamine 2,3-dioxygenase specific, cytotoxic T cells as immune regulators. Blood, 117(7): 2200-2210 (2011)

 **For updated literature or any other information please contact your local supplier.**

### Symbols:

	Storage temperature		Manufacturer		Contains sufficient for <n> tests
	Expiry date	<b>LOT</b>	Batch code	<b>RUO</b>	For research use only!
	Consult instructions for use	<b>CONT</b>	Content	<b>CE</b>	CE labelled
	Caution	<b>REF</b>	Catalogue number		



## Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident. Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

*For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.*

distributed in the US/Canada by:



**EAGLE**  
BIOSCIENCES

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH  
03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)