



EAGLE  
BIOSCIENCES

# Mouse / Rat Adrenaline (Epinephrine) ELISA Assay Kit

Catalog Number:

ADU39-K01 (1 x 96 wells)

*For Research Use Only. Not for use in diagnostic procedures.*

*v. 1.0*

---

EAGLE BIOSCIENCES, INC.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

WWW.EAGLEBIO.COM



## 1. Introduction and Principle of the Test

[The Mouse / Rat Adrenaline \(Epinephrine\) ELISA Assay kit](#) is an Enzyme-linked immunosorbent assay used for the quantitative and very sensitive determination of adrenaline in biological samples including serum, plasma, tissue, and cell culture samples. The Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit is for research use only and should not be used in diagnostic procedures.

Catecholamine is the name of a group of aromatic amines (noradrenaline, adrenaline, dopamine, and their derivatives) which act as hormones and neurotransmitter, respectively. Adrenaline and noradrenaline are formed from dopamine. They act on the cardiac musculature and the metabolism (adrenaline) as well as on the peripheral circulation (noradrenaline) and help the body to cope with acute and chronic stress.

An increased production of catecholamines can be found with tumours of the chromaffine system (pheochromocytoma, neuroblastoma, ganglioneuroma). An increased or decreased concentration of the catecholamines can also be found with hypertension, degenerative cardiac diseases, schizophrenia and manic-depressive psychosis.

The Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit provides materials for the quantitative measurement of adrenaline in low concentrated samples and for small sample volumes. Adrenaline is extracted using a cis-diol-specific affinity gel and acylated to N-acyladrenaline and then converted enzymatically into N-acylmetanephrine.

The competitive Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit uses the microtitre plate format. Adrenaline is bound to the solid phase of the microtiter plate. Acylated catecholamine from the sample and solid phase bound catecholamine compete for a fixed number of antiserum binding sites. When the system is in equilibrium, free antigen and free antigen-antiserum complexes are removed by washing. The antibody bound to the solid phase catecholamine is detected by anti-rabbit IgG / peroxidase. The substrate TMB / peroxidase reaction is monitored at 450 nm. The amount of antibody bound to the solid phase catecholamine is inversely proportional to the catecholamine concentration of the sample.

## 2. Precautions

- **For research use only.**
- Some reagents contain sodium azide as preservative. Avoid skin contact.
- Material of animal origin used in the preparation of the kit has been obtained from animals certified as healthy but these materials should be handled as potentially infectious.
- Some components of this kit are containing hazardous reagents. These components are marked with the adequate hazard label.



### 3. Storage and Stability

- On arrival, store the Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit at 2 - 8 °C. Once opened the Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit is stable until its expiry date. For stability of prepared reagents refer to Preparation of Reagents.
- Do not use components beyond the expiration date shown on the Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit labels.
- Do not mix various lots of any kit component within an individual assay.

### 4. Contents of the Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit

#### Reagents for Sample Preparation:

- 4.1 **Extraction Plate** **EX-PLATE** 2 plates  
48 wells  
coated with boronate affinity gel
- 4.2 **Extraction-Buffer** **EX-BUFF** 2 vials  
6 ml, ready for use
- 4.3 **HCl** **HCL** 1 vial  
21 ml, ready for use  
0.025 M HCl
- 4.4 **Standards (A - F)** **CAL A - F** 6 vials  
Each 4 ml, ready for use  
Concentrations:

Standard	A	B	C	D	E	F
<b>Adrenaline</b> (ng/ml)	0	0.15	0.5	1.5	5	25
<b>Adrenaline</b> (nmol/l)	0	0.82	2.7	8.2	27.3	137

- 4.5 **Control 1 & 2** **CON 1 & 2** 2 vials  
Each 4 ml, ready for use  
Concentrations: see q.c. certificate
- 4.6 **Acylation Reagent** **ACYL-REAG** 1 vial  
6 ml, ready for use  
Contains DMSO and DMF  
(please note that solvent reacts with many plastic materials including plastic trays; it does not react with normal pipette tips and with glass devices).
- 4.7 **Acylation Buffer** **ACYL-BUFF** 1 vial  
20 ml, ready for use



4.8	<b>Enzyme</b> each 2 ml, lyophilized Catechol-O-methyltransferase	<b>ENZYME</b>	3 vials
4.9	<b>Coenzyme</b> 1 ml, ready for use S-adenosyl-L-methionine	<b>COENZYME</b>	1 vial
4.10	<b>Enzyme Buffer</b> 3.5 ml, ready for use	<b>ENZYME-BUFF</b>	1 vial
4.11	<b>Enzyme Plate</b> 96 wells, ready for use	<b>ENZYME-PLATE</b>	1 piece
4.12	<b>Sample Stabilizer</b> 20 ml, ready for use	<b>STABILIZER</b>	1 vial

**Reagents for ELISA:**

4.13	<b>Adrenaline-Antiserum</b> 2.5 ml, ready for use, rabbit colour coded blue	<b>AS-AD</b>	1 vial
4.14	<b>MT-Strips</b> 8 wells each, break apart, precoated with adrenaline, colour coded blue	<b>STRIPS-AD</b>	12 strips
4.15	<b>POD Conjugate</b> 12 ml, ready for use, Anti-rabbit IgG-POD conjugate / peroxidase	<b>CONJ</b>	1 vial
4.16	<b>Wash Buffer</b> 20 ml, concentrate Dilute content with dist. water to 500 ml total volume	<b>WASH</b>	2 vials
4.17	<b>Substrate</b> 12 ml TMB solution, ready for use	<b>SUB</b>	1 vial
4.18	<b>Stop Solution</b> 12 ml, ready for use Contains 0.3 M sulphuric acid	<b>STOP</b>	1 vial
4.19	<b>Adhesive Foil</b> Ready for use	<b>FOIL</b>	10 pieces



Additional materials and equipment required but not provided:

- Pipettes (20, 50, 100, 150, 175, 280  $\mu$ l)
- Repeating dispenser for 20, 25, 50, 100, 150 and 1 ml
- Orbital shaker
- Multichannel pipette or Microplate washing device
- Microplate photometer
- Distilled water
- Heating cabinet with 37 °C (optional)

## 5. Sample Collection and Storage

### Plasma

- EDTA plasma samples are required for the assay. Physical and psychical stress usually causes a high increase of the catecholamine concentration. Therefore, it is recommended to let the patient rest for 20 to 30 minutes after the venipuncture and before collecting the blood sample.
- Haemolytic and especially lipemic samples should not be used for the assay, because false low values will be obtained with such samples.
- Immediately after collection the plasma samples should be centrifuged (preferable at 2 - 8 °C) and frozen. The samples are stable up to 1 week at -20 °C.
- To improve the stability each sample should be enriched with the Sample Stabilizer **STABILIZER** before freezing (max. 20% of the sample volume), e.g.:

Sample volume	Stabilizer volume
20 $\mu$ l	4 $\mu$ l
50 $\mu$ l	10 $\mu$ l
100 $\mu$ l	20 $\mu$ l
200 $\mu$ l	40 $\mu$ l
300 $\mu$ l	60 $\mu$ l
500 $\mu$ l	100 $\mu$ l

### Cell culture samples and various biological samples:

- The stability of such samples depends on the sample type and the way of collection. Therefore, a general procedure for collection and storage is not possible. However, it is recommended to freeze the samples immediately after collection. The samples should be stable at -20 °C for up to 1 week.



- To improve the stability each sample should be enriched with the Sample Stabilizer **STABILIZER** before freezing (max. 10% of the sample volume), e.g.:

Sample volume	Stabilizer volume
20 µl	2 µl
50 µl	5 µl
100 µl	10 µl
200 µl	20 µl
300 µl	30 µl
500 µl	50 µl

Acidified samples, which have a pH value of 5 or less must not be enriched with the Sample Stabilizer and have to be frozen immediately after collection.

### Tissue samples

Tissue samples can be homogenized in 0.01 N HCl in the presence of 0.15 mM EDTA and 4 mM sodium metabisulfite.

The following basic principles should be followed:

- Avoid excess of acid. This might exceed the buffer capacity of the extraction buffer. After adding the extraction buffer a pH value of 7 or above is mandatory. If the pH value is below 7 it is necessary to repeatedly add 50 µl of Extraction Buffer until the pH value is at or above 7.  
Acidified samples, which have a pH value of 5 or less must not be enriched with the Sample Stabilizer.
- Avoid substances in the samples with a cis-diol-structure (boric acid, Sorbitol, mannitol, etc.). These substances reduce the recovery of extraction.

## 6. Preparation of Reagents and Samples

### 6.1. Preparation of Reagents

#### 6.1.1 Wash Buffer

Dilute the contents of the bottle with distilled water to a total volume of 500 ml. For further use the diluted wash buffer must be stored at 2 - 8 °C for a maximum period of 4 weeks.

#### 6.1.2 Enzyme Mix

**NOTE:** The enzyme mix has to be prepared freshly prior to the assay (not longer than 10 - 15 minutes in advance). After use the reagent has to be discarded.

Reconstitute the content of one vial labelled **ENZYME** with 2 ml distilled water. Add 0.3 ml **COENZYME** and 0.7 ml **ENZYME-BUFF**, mix thoroughly (total volume: 3 ml).



The two additional bottles of **ENZYME** are allowing a second and a third run of the test. If the whole kit is to be used in one run it is sufficient to prepare one vial of enzyme mix.

All other reagents are ready for use.

## 6.2 Preparation of Samples

Allow reagents to reach room temperature.

Determinations in duplicates are recommended.

Each 20  $\mu$ l of Standards and Control 1 & 2 are extracted.

Each 1  $\mu$ l - 300  $\mu$ l of samples are extracted (alternatively: > 300  $\mu$ l up to 500  $\mu$ l).

1. Pipette each 20  $\mu$ l Standard A - F, 20  $\mu$ l Control 1 & 2 and each 1  $\mu$ l - 300  $\mu$ l Sample into the respective wells of the extraction plate.  
Correction for volume:  
Pipette 280  $\mu$ l of distilled water into the wells of the standards and controls (final volume: 300  $\mu$ l).  
Pipette as much distilled water into the wells of the samples to obtain a final volume of 300  $\mu$ l, e.g. 100  $\mu$ l sample + 200  $\mu$ l distilled water.  
For sample volumes above 300  $\mu$ l up to 500  $\mu$ l: fill up all wells to a final volume of 500  $\mu$ l.  
Within a run the final volume has to be the same in all wells (300  $\mu$ l or 500  $\mu$ l, respectively).
2. Pipette each 100  $\mu$ l Extraction Buffer into all wells.
3. Cover the plate with adhesive foil and incubate for 60 minutes at room temperature on an orbital shaker (high shaking rate).
4. Decant the plate and remove residual liquid by tapping the inverted plate on a paper towel.
5. Pipette each 1 ml prepared Wash Buffer into all wells and incubate for 5 minutes at room temperature on an orbital shaker (slow shaking rate).
6. Decant the plate and remove residual liquid by tapping the inverted plate on a paper towel.
7. Pipette each 150  $\mu$ l Acylation Buffer into all wells.
8. Pipette each 50  $\mu$ l Acylation Reagent into all wells and continue with step 9. immediately.  
(please note that solvent reacts with many plastic materials including plastic trays; it does not react with normal pipette tips and with glass devices)
9. Incubate for 20 minutes at room temperature on an orbital shaker (medium shaking rate).
10. Decant the plate and remove residual liquid by tapping the inverted plate on a paper towel.
11. Pipette each 1 ml prepared Wash Buffer into all wells and incubate for 5 minutes at room temperature on an orbital shaker (slow shaking rate).



12. Decant the plate and remove residual liquid by tapping the inverted plate on a paper towel.
13. Repeat the wash steps 11. and 12.
14. Pipette each 125  $\mu$ l HCl (0.025 M) for elution into all wells.
15. Cover the plate with adhesive foil and incubate for 20 minutes at room temperature on an orbital shaker (medium shaking rate).  
Caution: Do not decant the supernatant thereafter.
16. Transfer 100  $\mu$ l from the extraction plate into the respective wells of the enzyme plate.
17. Pipette each 20  $\mu$ l of freshly prepared Enzyme Mix (s. 6.1.2) into all wells of the enzyme plate. Colour changes to red.
18. Cover the plate with adhesive foil and incubate for 1 minute at room temperature on an orbital shaker (medium shaking rate).
19. Incubate the plate for 90 minutes at 37°C without shaking.  
(Alternatively: 120 minutes at room temperature (20 - 25°C) on an orbital shaker at medium shaking rate).  
Caution: Do not decant the supernatant thereafter.

Take each 100  $\mu$ l of the supernatant for the Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit.

## 7. Test Procedure ELISA

1. Pipette each 100  $\mu$ l prepared Standards, Controls and Samples into the respective wells (colour coded blue).
2. Pipette each 20  $\mu$ l Adrenaline-Antiserum (colour coded blue) into all wells.
3. Cover the plate with adhesive foil, shake briefly and incubate for 15 – 20 hours (overnight) at 2 - 6 °C.
4. Discard or aspirate the contents of the wells and wash thoroughly with each 250  $\mu$ l prepared Wash Buffer. Remove residual liquid by tapping the inverted plate on clean absorbent paper. Repeat the washing procedure 3 times.
5. Pipette each 100  $\mu$ l POD-Conjugate into all wells.
6. Incubate for 60 minutes at room temperature on an orbital shaker (medium shaking rate).
7. Washing: Repeat wash step 4.
8. Pipette each 100  $\mu$ l Substrate into all wells.





9. Incubate for 35 to 45 minutes at room temperature (20 – 25 °C) on an orbital shaker (medium shaking rate). Avoid exposure to direct sun light.
10. Pipette 100 µl Stop Solution into all wells.
11. Read the optical density at 450 nm (reference wavelength between 570 and 650 nm) in a microplate photometer within 15 minutes.

### 8. Calculation of Results

On a semilogarithmic graph paper the concentration of the standards (x-axis, logarithmic) are plotted against their corresponding optical density (y-axis, linear). Alternatively, the optical density of each standard and sample can be related to the optical density of the zero standard, expressed as the ratio OD/OD<sub>max</sub>, and then plotted on the y-axis.

A good fit is provided with 4 Parameter Logistic (alternatively Log-Logit or Cubic Spline).

The concentration of the controls can be read off the standard curve directly without any further conversion. The read concentrations of the samples have to be divided by a correction factor due to the use of 1 µl - 300 µl sample volume in relation to 20 µl standard.

$$\text{Correction factor} = \frac{\text{Sample volume for extraction } (\mu\text{l})}{20 \mu\text{l (Standard volume)}}$$

Example:  
300 µl sample was extracted and the concentration read off from the standard curve is 0.6 ng/ml.

$$\text{Correction factor} = 300 \mu\text{l} / 20 \mu\text{l} = 15$$

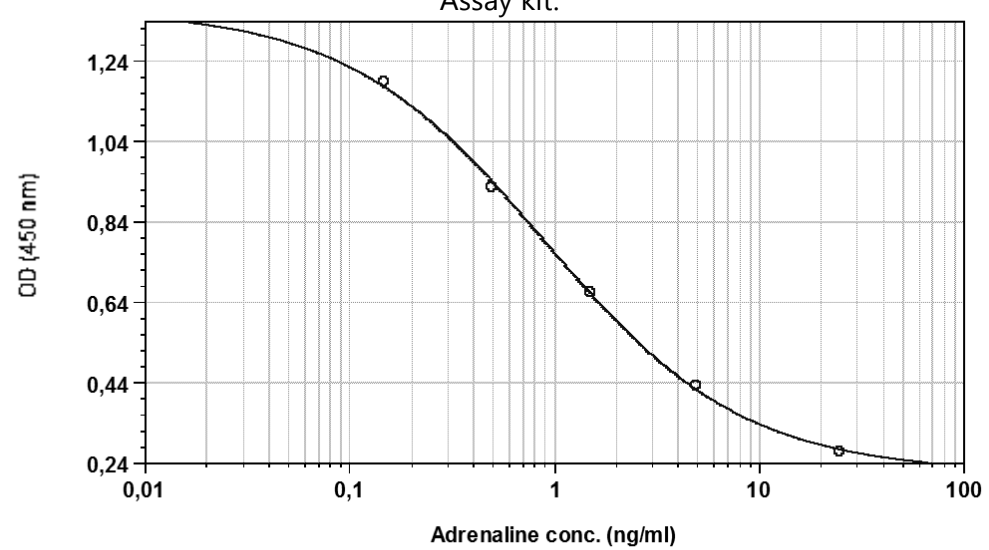
$$\text{Concentration of the sample} = 0.6 \text{ ng/ml} / 15 = 0.040 \text{ ng/ml} = 40 \text{ pg/ml}$$

Conversion into pmol/l:

$$\text{Adrenaline: } 1 \text{ pg} / \text{ml} = 5,46 \text{ pmol} / \text{l}$$

### Typical Example

Below a typical example of a standard curve with Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit:





## 9. Assay Characteristics

### 9.1 Reference Ranges

The reference ranges given below should only be taken as a guideline. It is recommended that each laboratory should establish its own normal values.

Plasma	< 100 pg/ml
--------	-------------

### 9.2 Sensitivity

The lower limit of detection of this Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit was determined by taking the 2fold standard deviation of the absorbance of the Zero Reference and reading the corresponding value from the standard curve. The sensitivity depends on the sample volume and can be calculated with the corresponding correction factor ( see 8. Calculation of Results)

Sensitivity:	$\frac{24 \text{ pg/ml (131 pmol/l)}}{\text{Correction factor}}$
Example for 300 $\mu$ l Sample (Correction factor 15):	$\frac{24 \text{ pg/ml}}{15} = 1.6 \text{ pg/ml}$ (8.7 pmol/l)

### 9.3 Specificity (Cross Reactivity)

Structural related components were tested for possible interference:

Substance	Cross Reactivity (%) Noradrenaline-Ab
Adrenaline	100
Noradrenaline	0.030
Dopamine	<0.01
Metanephrine	0.48
Normetanephrine	<0.01
3-Methoxytyramine	< 0.01
L-Dopa	< 0.002
Tyramine	< 0.002
Tyrosine	< 0.002
Homovanillic acid	< 0.002
Vanilmandelic acid	< 0.002



## 9.4 Recovery

Increasing amounts of adrenaline were added to an EDTA plasma sample and to a cell culture medium (RPMI 1640). Each spiked sample was assayed with the Mouse / Rat Adrenaline (Epinephrine) ELISA Assay kit. The analytical recovery was estimated at different concentrations by using the theoretically expected and the actually measured values.

Concentrations in pg/ml

EDTA-Plasma				cell culture medium			
added	measured	expected	% recovery	added	measured	expected	% recovery
0.0	4.5			0.0	5.1		
8.8	16.6	13.4	<b>124</b>	8.8	14.3	14.0	<b>103</b>
15.2	18.1	19.7	<b>92</b>	15.2	25.6	20.3	<b>126</b>
22.4	24.9	26.9	<b>93</b>	22.4	29.2	27.5	<b>106</b>
34.4	34.8	38.9	<b>89</b>	34.4	42.1	39.5	<b>107</b>
45.5	43.9	50.0	<b>88</b>	45.5	50.8	50.6	<b>100</b>
88.2	83.5	92.8	<b>90</b>	88.2	90.1	93.4	<b>97</b>
151.5	127.9	156.0	<b>82</b>	151.5	148.9	156.6	<b>95</b>
223.9	219.8	228.4	<b>96</b>	223.9	229.2	229.0	<b>100</b>
384.6	333.1	389.1	<b>86</b>	384.6	355.7	389.7	<b>91</b>
		mean:	<b>93</b>			mean:	<b>103</b>

## 9.5 Reproducibility

### Intra-Assay

The reproducibility of the ELISA method was investigated by determining the intra-assay-coefficients of variation (cv) by repeated measurements for EDTA-Plasma and cell culture medium (RPMI 1640) with different concentrations.

Concentrations in pg/ml

sample	n	mean value	sd	cv (%)
EDTA-Plasma	16	116.4	7.12	<b>6.1</b>
Cell culture medium	24	34.6	3.16	<b>9.2</b>



### Pipetting Scheme Sample Preparation

		Standards	Controls	Samples
Standard A - F	μl	20		
Control 1 & 2	μl		20	
Sample	μl			1 - 300
Dist. Water	μl	280	280	fill up to 300
Extraction Buffer	μl	100	100	100

Cover the plate with adhesive foil; shake for 60 minutes at room temperature  
Decant plate and remove residual liquid

Wash Buffer	ml	1	1	1
-------------	----	---	---	---

Shake for 5 minutes at room temperature (slow shaking rate)  
Decant plate and remove residual liquid

Acylation Buffer	μl	150	150	150
Acylation Reagent	μl	50	50	50

**Immediately:** Shake for 20 minutes at room temperature  
Decant plate and remove residual liquid

Wash Buffer	ml	1	1	1
-------------	----	---	---	---

Shake for 5 minutes at room temperature (slow shaking rate)  
Decant plate and remove residual liquid

Wash Buffer	ml	1	1	1
-------------	----	---	---	---

Shake for 5 minutes at room temperature (slow shaking rate)  
Decant plate and remove residual liquid

HCl	μl	125	125	125
-----	----	-----	-----	-----

Cover the plate with adhesive foil; Shake for 20 min at room temperature  
**Caution: Do not decant the supernatant thereafter**

Transfer Enzyme Plate	to μl	100	100	100
-----------------------	-------	-----	-----	-----

Enzyme Mix (fresh)	μl	20	20	20
--------------------	----	----	----	----

Cover the plate with adhesive foil; Shake for 1 minute at room temperature  
Incubate for 90 minutes at 37°C

**Caution: Do not decant the supernatant thereafter**

For the ELISA transfer each 100 μl



### Pipetting Scheme ELISA

	Standards	Controls	Samples
--	-----------	----------	---------

<b>Adrenaline (blue)</b>				
Standard A - F	μl	100		
Controls 1 & 2	μl		100	
Samples	μl			100
Adrenaline Antiserum	μl	20	20	20

Cover the plates with adhesive foil and shake briefly  
Incubate for 15 – 20 hours (overnight) at 2 - 6 °C

4 x washing

POD-Conjugate	μl	100	100	100
---------------	----	-----	-----	-----

Shake for 60 minutes at room temperature

4 x washing

Substrate	μl	100	100	100
-----------	----	-----	-----	-----

Shake for 35 – 45 minutes at room temperature

Stop Solution	μl	100	100	100
---------------	----	-----	-----	-----

Reading of absorbance at 450 nm



## 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, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.*