



## **Rat Tissue-type Plasminogen Activator (tPA) Active ELISA**

Catalog Number: TPA29-K01  
96 Wells  
For Research Use Only  
*v. 1.0*

---

EAGLE BIOSCIENCES, INC.  
20A NW Blvd., Suite 112, Nashua, NH 03063  
Phone: 866-419-2019 Fax: 617-419-1110  
[WWW.EAGLEBIO.COM](http://WWW.EAGLEBIO.COM)



---

## INTRODUCTION

---

Tissue-type Plasminogen Activator (tPA) is a member of the serine proteinase family. tPA functions to lyse fibrin clots into soluble plasmin fragments. tPA is active in two forms, single-chain and two-chain. The two-chain tPA is created via interaction with the plasmin product cleaving the single chain. This two-chain form is regarded as the more active form.

Both single-chain and two-chain tPA are complexable with PAI-1. PAI-1 acts as an inhibitor for tPA by binding to the tPA and thus stifling its ability to lyse fibrin. tPA can serve as an indicator of both myocardial infarction for patients with impaired fibrinolytic systems as well as a marker for type-II diabetes.

---

## PRINCIPLES OF PROCEDURE

---

This Rat Tissue-type Plasminogen Activator (tPA) Active ELISA (Enzyme-Linked Immunosorbent Assay) is for the quantitative analysis of active tPA levels in biological fluid. This test kit operates on the basis of sandwich ELISA where free, active, tPA enzyme complexes with PAI-1 and is quantified with the use of an HRP labeled secondary antibody.

First the biotinylated PAI-1 binds to the avidin coated wells. Next, active tPA present in the standard or unknown, complexes with PAI-1. Inactive or complexed tPA is removed in a subsequent wash step. A primary antibody specific for tPA is then added to each well followed by the HRP conjugated secondary antibody. The bound conjugated secondary antibody is detected by the addition of substrate, which generates an optimal color after 10 minutes. Quantitative test results may be obtained by the measure and comparison of the sample and standard absorbance readings when read with a microplate reader at 450 nm.

---

## MATERIALS PROVIDED

---

Component	Contents	Quantity	Storage
<b>Coated Plate</b>	Coated 96-well plate	1 plate	4°C
<b>Rat tPA Activity Standard</b>	Lyophilized active rat tPA standard	1 vial	4°C
<b>Biotinylated PAI-1</b>	Lyophilized biotinylated PAI-1	1 vial	4°C
<b>10x Wash Buffer</b>	10x Solution for washing plate	50 mL	4°C
<b>Substrate</b>	TMB substrate	10 mL	4°C
<b>Primary Antibody</b>	Lyophilized anti-mouse tPA antibody	1 vial	4°C
<b>Secondary Antibody</b>	HRP-conjugated secondary antibody	1 vial	4°C

---

## MATERIALS NEEDED BUT NOT PROVIDED

---

1. Adjustable pipettes (10-1,000 µL) and disposable tips
2. Beakers, flasks, and cylinders as necessary for preparation of reagents
3. Microplate reader with 450 nm filter
4. Deionized Water
5. 1 N Sulfuric Acid
6. TBS Buffer (see Reagent Preparation)
7. Blocking Buffer (see Reagent Preparation)
8. Microplate shaker with uniform horizontal circular movement up to 300 rpm
9. Plastic film or plate cover to cover plate during incubation



---

## STORAGE CONDITIONS

---

1. Store this kit and its components at 4°C until use.
2. Do not freeze.

---

## WARNINGS AND PRECAUTIONS

---

1. Use aseptic technique when opening and dispensing reagents.
2. This kit is designed to work properly as provided and instructed. Additions, deletions, or substitutions to the procedure or reagents are not recommended, as they may be detrimental to the assay.
3. Exercise universal precautions during the performance or handling of this kit or any component contained therein.

---

## PROCEDURAL NOTES

---

1. This assay should be performed at room temperature.
2. Performance of the entire kit at once is not required. When performing this kit in part, please adhere to the following:
  - All unused components should be returned to storage at 4°C.
  - Desiccant bag must remain in foil pouch with unused strips. Keep pouch sealed when not in use to maintain a dry environment. Seal with a heat sealer. If a heat sealer is not available, thoroughly close the open end with tape. Try to remove excess air before sealing.
  - If not using the entire plate at once, then prepare only the required amount of reagents where appropriate.
  - tPA Standard and reconstituted biotinylated PAI-1 stock solutions must be stored at -70 °C after initial use.
3. To minimize error due to handling, wipe the exterior bottom of the microplate wells with a lint-free paper towel.

---

## SAMPLE COLLECTION AND PREPARATION

---

Samples should be collected using trisodium citrate, acidified citrate or Stabilite™ (DiaPharma) collection media. Collection should be in accordance with the collection vials manufacturers instructions or in a 1:10 ratio of collection media to blood.

Immediately, upon collection of blood, the samples should be centrifuged at 3000 x g. This should ensure the removal of platelets as they can release PAI-1 that in turn complexes with uPA. The plasma can be transferred to a clean plastic tube and stored frozen for up to one month.

Samples are stable for approximately 5 hours when stored at 4° C with the Sabilyte™ media or up to one month when stored at -20°C.

**Note:** Detergents such as Triton X cause interference with the assay. If using detergent extracted samples, it is necessary to dialyze the samples overnight to remove the detergent.

---

## REAGENT PREPARATION

---

The following solutions should be prepared fresh before starting the assay.

1. **TBS Buffer:** 0.10 M TRIS, 0.15 M NaCl, pH 7.4
2. **Blocking Buffer:** 3% BSA in TBS Buffer.
3. **Wash Buffer:** Dilute the 10x Wash Buffer to 1x with deionized water and mix prior to use.



---

## ASSAY PROCEDURE

---

1. Remove microplate from the bag.
2. Make a working solution of Biotinylated PAI-1 by reconstituting according to the amount indicated on the vial using 3% BSA Blocking Buffer.
3. Add 100  $\mu$ L of the Biotinylated PAI-1 Solution to both the standard and test wells.
4. Shake the plate at 300 rpm on a plate shaker for 30 minutes.
5. Wash wells according to the following wash procedure:
  - a. Remove contents of the plate by inversion into an appropriate disposal device.
  - b. Tap out remaining contents of the plate onto a lint free paper towel.
  - c. Add 300  $\mu$ L of Wash Buffer.
  - d. Let stand for 2-3 minutes.
  - e. Repeat procedure 2 more times and then proceed to step “f”.
  - f. Remove contents of the plate by inversion into an appropriate disposal device.
  - g. Tap out the remaining contents of the plate onto a lint free paper towel then proceed to step 6.

**Note:** The decanted wells should be void of visible moisture before proceeding. If moisture is still visible then follow step “g” until satisfactory results are obtained.

6. Prepare the Standards as indicated in the provided dilution table.

**Note:** The standards should be applied to the plate immediately upon preparation.
7. Add 100  $\mu$ L of Standards or unknowns to the plate. See **Scheme I** for suggested template design.
8. Shake plate at 300 rpm on the plate shaker for 30 minutes.
9. Wash wells according to step 5 located above in this section.
10. Make a working solution of Primary Antibody by reconstituting according to the amount indicated on the vial using 3% BSA Blocking Buffer.
11. Add 100  $\mu$ L of the Primary Antibody Solution to each well.
12. Shake plate at 300 rpm on the plate shaker for 30 minutes.
13. Wash wells according to step 5 located above in this section.
14. Make a working solution of Secondary Antibody by reconstituting according to the amount indicated on the vial using 3% BSA Blocking Buffer.
15. Add 100  $\mu$ L of the Secondary Antibody Solution to each well.
16. Shake plate at 300 rpm on the plate shaker for 30 minutes.
17. Wash wells according to step 5 located above in this section.
18. Add 100  $\mu$ L of TMB Substrate to each well and incubate for 10 minutes.
19. Stop the reaction with 50  $\mu$ L per well of 1 N  $H_2SO_4$  and read plate at 450 nm.
20. If accounting for substrate background, use 2 to 8 wells as blanks with only substrate in the wells (100  $\mu$ L/well). Subtract the average of these absorbance values from the absorbance values of the wells being assayed.

**NOTE:** Some microplate readers can be programmed to do these subtractions automatically when reading the plate. Consult your instrument manual.



### Scheme I: Suggested Plate Layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	S <sub>8</sub>	S <sub>7</sub>	S <sub>6</sub>	S <sub>5</sub>	S <sub>4</sub>	S <sub>3</sub>	S <sub>2</sub>	S <sub>1</sub>	B <sub>0</sub>	U <sub>1</sub>	U <sub>2</sub>	U <sub>3</sub>
B	S <sub>8</sub>	S <sub>7</sub>	S <sub>6</sub>	S <sub>5</sub>	S <sub>4</sub>	S <sub>3</sub>	S <sub>2</sub>	S <sub>1</sub>	B <sub>0</sub>	U <sub>1</sub>	U <sub>2</sub>	U <sub>3</sub>
C	U <sub>4</sub>	U <sub>5</sub>	U <sub>6</sub>	U <sub>7</sub>	U <sub>8</sub>	U <sub>9</sub>	U <sub>10</sub>	U <sub>11</sub>	U <sub>12</sub>	U <sub>13</sub>	U <sub>14</sub>	U <sub>15</sub>
D	U <sub>4</sub>	U <sub>5</sub>	U <sub>6</sub>	U <sub>7</sub>	U <sub>8</sub>	U <sub>9</sub>	U <sub>10</sub>	U <sub>11</sub>	U <sub>12</sub>	U <sub>13</sub>	U <sub>14</sub>	U <sub>15</sub>
E	U <sub>16</sub>	U <sub>17</sub>	U <sub>18</sub>	U <sub>19</sub>	U <sub>20</sub>	U <sub>21</sub>	U <sub>22</sub>	U <sub>23</sub>	U <sub>24</sub>	U <sub>25</sub>	U <sub>26</sub>	U <sub>27</sub>
F	U <sub>16</sub>	U <sub>17</sub>	U <sub>18</sub>	U <sub>19</sub>	U <sub>20</sub>	U <sub>21</sub>	U <sub>22</sub>	U <sub>23</sub>	U <sub>24</sub>	U <sub>25</sub>	U <sub>26</sub>	U <sub>27</sub>
G	U <sub>28</sub>	U <sub>29</sub>	U <sub>30</sub>	U <sub>31</sub>	U <sub>32</sub>	U <sub>33</sub>	U <sub>34</sub>	U <sub>35</sub>	U <sub>36</sub>	U <sub>37</sub>	U <sub>38</sub>	U <sub>39</sub>
H	U <sub>28</sub>	U <sub>29</sub>	U <sub>30</sub>	U <sub>31</sub>	U <sub>32</sub>	U <sub>33</sub>	U <sub>34</sub>	U <sub>35</sub>	U <sub>36</sub>	U <sub>37</sub>	U <sub>38</sub>	U <sub>39</sub>

### CALCULATIONS

1. Subtract the average O.D. value of the blank wells (BLK) from all other pairs of wells.
2. Average the O.D. values for each pair of duplicate wells.
3. Plot a standard curve using the average O.D. value for each standard value versus the concentration of standard.
4. Determine the concentration of each unknown by interpolation from the standard curve.

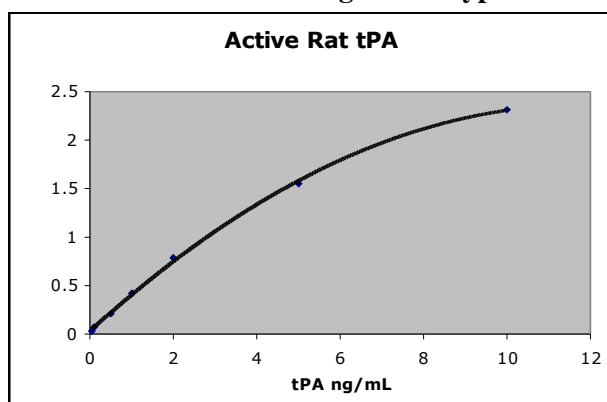
### TYPICAL RESULTS

**Note:** "Typical data" is a representation. Variances in data will occur. Optical density readings may fluctuate during the shelf-life or due to lot variance.

#### Typical Data:

Standard	[tPA] (ng/mL)	O. D.
S <sub>1</sub>	0.05	0.030
S <sub>2</sub>	0.1	0.030
S <sub>3</sub>	0.25	0.071
S <sub>4</sub>	0.5	0.212
S <sub>5</sub>	1.0	0.423
S <sub>6</sub>	2.0	0.789
S <sub>7</sub>	5.0	1.553
S <sub>8</sub>	10.0	2.310

**Figure 1: Typical Standard Curve**





---

## PERFORMANCE CHARACTERISTICS

---

Assay Range: 0.05- 10 ng/mL

Samples with tPA levels higher than 10 ng/mL should be diluted in similar media devoid of active tPA.

---

## REFERENCES

---

1. Thogersen A et al. (1998) Circulation 98: 2241-2247
2. Eliasson M et al. (2003) Cardiovascular Diabetology 2:19

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

*Product Developed and Manufactured in the USA*