



Plasminogen (human) ELISA kit

Catalog # ADI-900-231
96-Well Kit



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Please contact Enzo Life Sciences Technical Support if necessary.



Carefully note the handling and storage conditions of each kit component.

TABLE OF CONTENTS

Introduction	2
Principles of Method	2
Intended Use	2
Storage and Stability	3
Chemical Hazard	3
Kit Contents	3
Other Materials Needed	4
Reagent Preparation	5
Human Plasminogen standard	5
Secondary Antibody	5
AV-HRP	6
Washing buffer	6
Sample preparation	7
Assay Procedure	7
Characteristics	9
Typical result	9
Sensitivity	10
Specificity	10
Precision	10
Recovery	11
Troubleshooting	12
References	13
Contact Information	14

INTRODUCTION

Plasminogen, a 92kDa glycoprotein, is produced by the liver and is present in plasma and extracellular fluids. Plasminogen is the inactive precursor of plasmin, a potent serine protease involved in the dissolution of fibrin blood clots. Plasminogen can be converted into the active plasmin by plasminogen activators urokinase (uPA), tissue plasminogen activator (tPA), factor XII-dependent components. The plasmin system has been implicated in a variety of physiological and pathological processes such as fibrinolysis, tissue remodeling, cell migration, inflammation, and tumor invasion and metastasis. Hereditary defects of plasminogen are a predisposing risk factor for thromboembolic disease.

PRINCIPLES OF METHOD

The design of this assay is based on a sandwich Enzyme-Linked Immunosorbent Assay (ELISA). The microtiter plate provided in this kit has been pre-coated with a monoclonal antibody specific to human plasminogen. Samples are pipetted into these wells. Nonbound plasminogen and other components of the sample should be removed by washing, then biotin-conjugated monoclonal antibody specific to plasminogen added. In order to quantitatively determine the amount of plasminogen present in the sample, Avidin conjugated to Horseradish Peroxidase (HRP) should be added to each microplate well. The final step, a TMB-substrate solution added to each well. Finally, a sulfuric acid solution is added and the resulting yellow colored product is measured at 450nm. Since the increases in absorbency is directly proportional to the amount of captured plasminogen.

INTENDED USE

The Enzo Life Sciences Plasminogen (human) ELISA kit is to be used for the *in vitro* quantitative determination of human plasminogen in human serum, human plasma, cell lysate and buffered solution. The assay will recognize native human plasminogen.

This kit has been configured for research use only and is not to be used in diagnostic procedures.

STORAGE AND STABILITY

All kit components of this kit are stable at 2-8°C. Any unused reconstituted standard should be discarded or frozen at -70°C. **Standard can be frozen and thawed one time only without loss of immunoreactivity.**

CHEMICAL HAZARD

Stop solution: This reagent is an irritant to eyes, skin and mucous membranes. Avoid contact with eyes, skin and clothing. Wear suitable protective clothing, gloves and eye protection. In the event of contact with eyes or skin, wash immediately with plenty of water.

Standard protein and 2nd Antibody containing Sodium Azide as a preservative.

KIT CONTENTS

Contents	Number	Volume
96 Well Plate	1 (in aluminum foil bag with desiccant)	
Incubation Buffer	1	30ml
Washing Buffer	2	(20X) 25ml
Standard Protein	1 Glass vial (lyophilized)	
Standard/Sample Dilution Buffer	1	25ml
Secondary Antibody	1 Glass vial (lyophilized)	
AV-HRP	1	150µl
Secondary Antibody/ AV-HRP Dilution Buffer	1	25ml
Substrate (TMB)	1	15ml
Stop Solution	1	15ml
Protocol booklet	1	-
Plate sealers	2	-

1. **96 Well Plate:**
Human plasminogen microtiter plate, one plate of 96 wells

A plate using break-apart strips coated with a mouse monoclonal antibody specific to human plasminogen.
2. **Standard Protein:**
Native human plasminogen.
3. **Secondary Antibody:**
Biotinylated anti human plasminogen.
4. **AV-HRP**
Avidin linked Horseradish Peroxidase (HRP, enzyme)
5. **Substrate** (Stabilized chromogen)
Tetramethylbenzidine (TMB) solution
6. **Stop Solution**
1N solution of sulfuric acid (H₂SO₄).
7. **Plate sealer**
Adhesive sheet.

Do not mix or interchange different reagents from various kit lots.

OTHER MATERIALS NEEDED

1. Microtiter plate reader capable of measurement at or near 450nm.
2. Calibrated, adjustable precision pipettes, preferably with disposable plastic tips (A manifold multi-channel pipette is desirable for large assays.)
3. Distilled or deionized water
4. Data analysis and graphing software
5. Vortex mixer
6. Polypropylene tubes for diluting and aliquoting standard
7. Absorbent paper towels
8. Calibrated beakers and graduated cylinders of various sizes

REAGENT PREPARATION

Human Plasminogen standard

1. Reconstitute the lyophilized Human Plasminogen standard by adding 1ml of **Standard/Sample Dilution Buffer** to make the 0.2 μ g/ml standard stock solution. Allow solution to sit at RT for 5 minutes, then gently vortex to mix completely. Use within one hour of reconstituting. Standard can be frozen and thawed one time only without loss of immunoreactivity.
2. Prepare 1ml of 64ng/ml top standard by adding 320 μ l of the above stock solution in 750 μ l of **Standard/Sample Dilution Buffer**. Perform 2-fold serial dilutions of the top standards to make the standard curve within the range of this assay (1ng/ml ~ 64ng/ml) as below. **Standard/Sample Dilution Buffer** serves as the zero standard (0ng/ml).

Standard	Add	Into
64ng/ml	320.0 μ l of the std. (0.2 μ g/ml)	680.0 μ l of the Standard/ Sample Dilution Buffer
32ng/ml	500 μ l of the std. (64ng/ml)	500.0 μ l of the Standard/ Sample Dilution Buffer
16ng/ml	500 μ l of the std. (32ng/ml)	500.0 μ l of the Standard/ Sample Dilution Buffer
8ng/ml	500 μ l of the std. (16ng/ml)	500.0 μ l of the Standard/ Sample Dilution Buffer
4ng/ml	500 μ l of the std. (8ng/ml)	500.0 μ l of the Standard/ Sample Dilution Buffer
2ng/ml	500 μ l of the std. (4ng/ml)	500.0 μ l of the Standard/ Sample Dilution Buffer
1ng/ml	500 μ l of the std. (2ng/ml)	500.0 μ l of the Standard/ Sample Dilution Buffer
0ng/ml	1.0ml of the Standard/Sample Dilution Buffer	

Secondary Antibody

100X secondary antibody solution can be made by adding 150 μ l secondary antibody/AV-HRP dilution buffer in the vial.

1. Equilibrate to room temperature, mix gently.
2. Mix 20 μ l **Secondary Antibody concentrated solution (100X)** + 1.98ml **Secondary Antibody/AV-HRP dilution buffer**.

Label as “*Working Secondary antibody Solution*”.

3. Return the unused **Secondary Antibody concentrated solution** to the refrigerator.

AV-HRP

1. Equilibrate to room temperature, mix gently.
2. Mix 20µl **AV-HRP concentrated solution** (100X) + 1.98ml **Secondary Antibody/AV-HRP dilution buffer**.
3. Label as “Working AV-HRP Solution”.
4. Return the unused **AV-HRP concentrated solution** to the refrigerator.

Washing buffer

1. Equilibrate to room temperature, mix to re-dissolve any precipitated salt.
2. Mix 0.5 volume **Wash buffer concentrate solution** (20X) + 9.5 volumes of deionized water. Label as “*Working Washing Solution*”.
3. Store both the concentrated and the Working Washing Solution in the refrigerator.

***Directions for washing**

1. Fill the wells with 300µl of “Working Washing Buffer”. Let soak for 1 to 3 minutes and then all residual wash-liquid must be drained from the wells by aspiration (taking care not to scratch the inside of the well) or decantation, followed by forceful tapping of the plate on absorbent paper. Never insert absorbent paper directly into the wells.

If using an automated washer, the operating instructions for washing equipment should be carefully followed.

2. Incomplete washing will adversely affect the assay and renders false results.
3. It is recommended to use laboratory tape to hold plate strips to the plate frame while performing the plate washing to avoid strips coming free of the frame.

SAMPLE PREPARATION

Blood should be collected by veinpuncture. For plasma samples, blood may be drawn into tubes containing sodium citrate or heparin, EDTA. The serum or plasma should be separated from the coagulated or packed cells by centrifugation. Specimens may be shipped at room temperature and then stored refrigerated at 2-8°C if testing is to take place within one week after collection. If testing is to take place later than one week, specimens should be stored at -20°C. Avoid repeated freeze/thawing. Serum and plasma require at least 2000 fold dilution in the Standard/Sample Dilution Buffer.

Assay Procedure

- Allow all reagents to reach room temperature before use. Gently mix all liquid reagents prior to use.
 - All standards, controls and samples should be run in duplicate for confirmation of reproducibility.
 - A standard curve must be run with each assay.
 - If particulate matter is present in the analyte, centrifuge or filter prior to analysis.
 - Maintain a consistent order of components and reagents addition from well to well. This ensures equal incubation times for all wells.
1. Determine the number of strips needed for assay. Insert these in the frame(s) for current use. (Re-bag extra strips and frame. Refrigerate for further use.)
 2. Add 300µl of Incubation buffer to all wells and incubate the plate for 5 minutes at room temperature.
 3. Thoroughly aspirate or decant the solution from the wells. Wash wells 2 times (See “*Directions for washing*”).
 4. For the standard curve, add 100µl of the standard to the appropriate microtiter wells. Add 100µl of the **Standard/Sample Dilution Buffer** to zero wells.
 5. Serum and plasma require at least 2000 fold dilution in the **Standard/Sample Dilution Buffer**. And add 100µl of sample to each well.
 6. Cover the plate with the plate cover and incubate for 2 hours at room temperature.
 7. Thoroughly aspirate or decant the solution from the wells. Wash the wells 3 times (See “*Directions for washing*”).
 8. Pipette 100µl of “Working Secondary Antibody Solution” into each well.

9. Cover the plate with the plate cover and incubate for 1 hour at room temperature.
10. Thoroughly aspirate or decant the solution from the wells. Wash the wells 3 times (See “*Directions for washing*”).
11. Add 100µl “Working AV-HRP Solution” to each well.
12. Cover the plate with the plate cover and incubate for 30 minutes at room temperature.
13. Thoroughly aspirate or decant the solution from the wells. Wash the wells 3 times (See “*Directions for washing*”).
14. Add 100µl of **Substrate** to each well. The liquid in the wells should begin to turn blue.
15. Incubate the plate at room temperature.

- Do not cover the plate with aluminum foil, or color may develop.

The incubation time for chromogen substrate is often determined by the microtiter plate reader used. O.D. values should be monitored and the substrate reaction stopped before O.D. of the positive wells exceeds the limits of the instrument. O.D. values at 450nm can only be read after the Stop Solution has been added to each well.

- Because the **Substrate** is light sensitive, avoid prolonged exposure to light for the remaining Substrate solution.
 - Typically, reaction is stopped 5~10 minutes after treatment of Substrate, but this time can be adjusted as the user desires.
16. Add 100µl of **Stop Solution** to each well. The solution in the wells should change from blue to yellow.
 17. Read the absorbance of each well at 450nm. Read the plate within 20 minutes of adding the Stop Solution.
 18. Plot the absorbance of the standard against the standard concentration (Optimally, the background absorbance can be subtracted from all data points, including standards, unknowns and controls, prior to plotting.). Draw a smooth curve through these points to construct the standard curve.
 19. Read the human plasminogen concentrations for the unknown samples and controls from the standard curve plotted in step 18. Multiply value(s) obtained for the unknown sample by the dilution factor (Samples producing signals greater than that of the highest standard should be further diluted in the **Standard/Sample Dilution Buffer**).

CHARACTERISTICS

Typical results

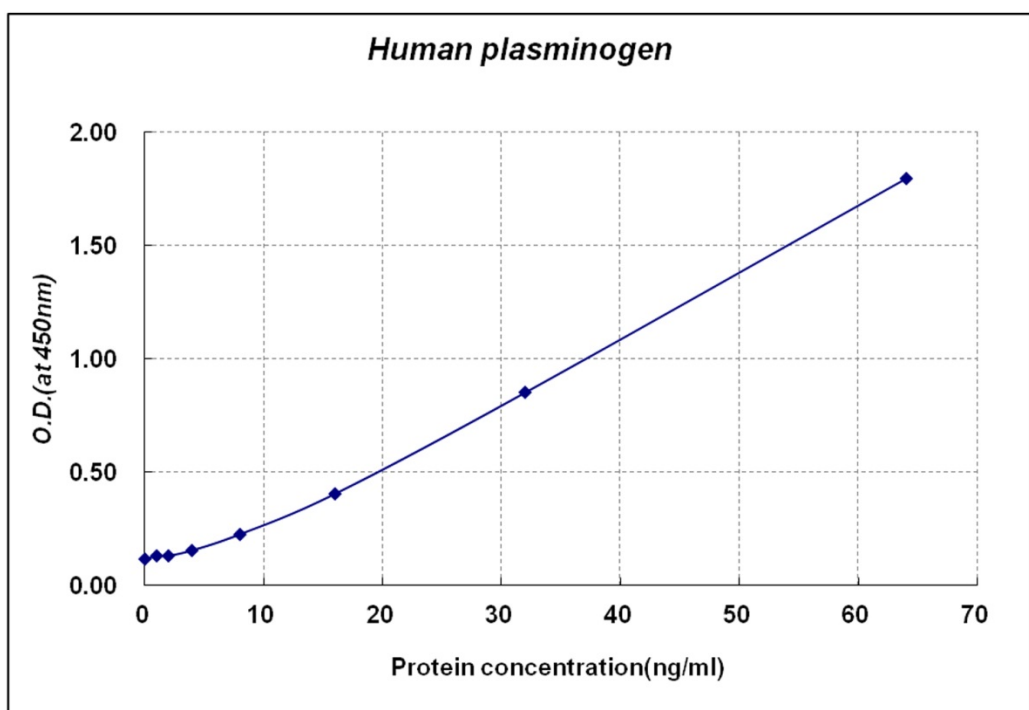
The standard curve below is for illustration only and should not be used to calculate results in your assay.

A standard curve must be run with each assay.

Standard Human Plasminogen (ng/ml)	Optical Density (at 450nm)
0	0.116
1	0.129
2	0.131
4	0.155
8	0.226
16	0.405
32	0.853
64	1.798

Limitations

- Do not extrapolate the standard curve beyond the 64ng/ml standard point.
- Other buffers and matrices have not been investigated.
- The rate of degradation of native human plasminogen in various matrices has not been investigated.



Sensitivity

The minimal detectable dose of human plasminogen was calculated to be 2.01ng/ml, by subtracting two standard deviations from the mean of 12 zero standard replicates (ELISA buffer, S0) and intersecting this value with the standard curve obtained in the same calculation.

N	1	2	3	4	5	6	7	8	9	10	11	12
ZERO	0.083	0.067	0.069	0.068	0.069	0.069	0.083	0.067	0.069	0.068	0.069	0.069

Average	SD	LLD	LLD mean (ng/ml)
0.0708	0.0057	0.0823	2.0136

Specificity

The following substances were tested and found to have no cross-reactivity: human serum albumin, transferrin, IgG, alpha-fetoprotein (AFP), fibrinogen, Vitamin D binding protein (VDBP), Hemoglobin.

Precision

1. Within-Run (Intra-Assay)

(n=6)

Mean (ng/ml)	SD (ng/ml)	CV (%)
6.23	0.59	9.47
15.05	0.64	4.27
33.36	0.91	2.73
63.91	1.56	2.44

2. Between-Run (Inter-Assay)

(n=4)

Mean (ng/ml)	SD (ng/ml)	CV (%)
6.52	0.14	2.18
15.29	0.61	4.00
34.0	0.60	1.78
63.50	0.23	0.37

Recovery

Recovery on addition is 97.12~103.00% (mean 100.14%).

- Plasminogen in Serum A: 102.177µg/ml;
- Plasminogen in Serum B: 114.834µg/ml

Analyte added (ng/ml)	Added analyte+1/256000 diluted Serum A (ng/ml)	Recovery (%)	Added analyte+1/256000 diluted Serum B (ng/ml)	Recovery (%)
8	8.399	103.0	8.495	101.10
16	16.399	100.12	16.495	99.15
32	32.399	97.17	32.495	100.56
64	64.399	102.60	64.495	97.42

For the recovery assay, a known amount of the ELISA standard is spiked into serially-diluted normal human serum and run in the assay. Any endogenous levels are taken into account and then % recovery of the exogenously spike standard is determined.

TROUBLESHOOTING

Problem	Possible Cause	Solution
High signal and background in all wells	Insufficient washing	Increase number of washes Increase time of soaking between in wash
	Too much AV-HRP	Check dilution, titration
	Incubation time too long	Reduce incubation time
	Development time too long	Decrease the incubation time before the stop solution is added
No signal	Reagent added in incorrect order, or incorrectly prepared	Review protocol
	Standard has gone bad (If there is a signal in the sample wells)	Check the condition of stored standard
	Assay was conducted from an incorrect starting point	Reagents allows to come to 20~30°C before performing assay
Too much signal – whole plate turned uniformly blue	Insufficient washing – unbound AV-HRP remaining	Increase number of washes carefully
	Too much AV-HRP	Check dilution
	Plate sealer or reservoir reused, resulting in presence of residual AV-HRP	Use fresh plate sealer and reagent reservoir for each step

Problem	Possible Cause	Solution
Standard curve achieved but poor discrimination between point	Plate not developed long enough	Increase substrate solution incubation time
	Improper calculation of standard curve dilution	Check dilution, make new standard curve
No signal when a signal is expected, but standard curve looks fine	Sample matrix is masking detection	More diluted sample recommended
Samples are reading too high, but standard curve is fine	Samples contain protein levels above assay range	Dilute samples and run again
Edge effect	Uneven temperature around work surface	Avoid incubating plate in areas where environmental conditions vary. Use plate sealer.

REFERENCES

1. Castellino F.J. and Ploplis V.A., 2005, Thromb Haemost. 93:647-654
2. Wun T.C., 1988, Crit Rev Biotechnol. 8:131-148



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com