

ab108893 – Human Plasminogen (PLG) ELISA Kit

Instructions for Use

For the quantitative measurement of Human Plasminogen (PLG) in plasma, serum, urine, saliva, milk and cerebrospinal fluid

This product is for research use only and is not intended for diagnostic use.

Table of Contents

INTRODUCTION

1.	BACKGROUND	1
2.	ASSAY SUMMARY	2

GENERAL INFORMATION

3.	PRECAUTIONS	3
4.	STORAGE AND STABILITY	3
5.	MATERIALS SUPPLIED	3
6.	MATERIALS REQUIRED, NOT SUPPLIED	4
7.	LIMITATIONS	4
8.	TECHNICAL HINTS	5

ASSAY PREPARATION

9.	REAGENT PREPARATION	6
10.	STANDARD PREPARATIONS	9
11.	SAMPLE PREPARATION	12
12.	PLATE PREPARATION	14

ASSAY PROCEDURE

13.	ASSAY PROCEDURE	15
-----	-----------------	----

DATA ANALYSIS

14.	CALCULATIONS	17
15.	TYPICAL DATA	18
16.	TYPICAL SAMPLE VALUES	19
17.	ASSAY SPECIFICITY	21

RESOURCES

18.	TROUBLESHOOTING	21
19.	NOTES	24

1. BACKGROUND

Abcam's Plasminogen (PLG) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Plasminogen concentrations in Human plasma, serum, urine, saliva, milk and cerebrospinal fluid.

A Plasminogen specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Plasminogen specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Plasminogen captured in plate.

Plasminogen is a single chain glycoprotein zymogen that is synthesized in the liver and circulated in plasma with a molecular weight of 90 kDa. The N-terminal portion of the molecule is made up of five kringle domains that bind to fibrin. The native molecule has an amino-terminal glutamic acid, known as glu-plasminogen, but this can undergo proteolytic cleavage by plasmin to lys-plasminogen. The inactive proenzyme Plasminogen is converted to the active enzyme plasmin that ultimately digests fibrin. Tissue-type Plasminogen activator (tPA) or urokinase-type Plasminogen activator (uPA) catalyzes the activation of Plasminogen, while Plasminogen activator inhibitors (PAIs) inhibits the activation. The Plasminogen system plays a role in macrophage recruitment, arterial stenosis, atherosclerosis, aneurysm formation, skin and corneal wound healing, glomerulonephritis, and neovascularization.

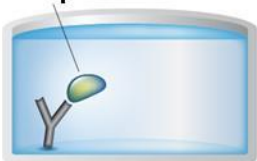
2. ASSAY SUMMARY

Primary capture antibody



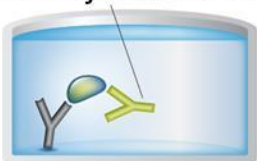
Prepare all reagents, samples and standards as instructed.

Sample



Add standard or sample to each well used. Incubate at room temperature.

Primary detector antibody



Wash and add prepared biotin antibody to each well. Incubate at room temperature.

Streptavidin Label



Wash and add prepared Streptavidin-Peroxidase Conjugate. Incubate at room temperature.

Substrate **Colored product**



Add Chromogen Substrate to each well. Incubate at room temperature. Add Stop Solution to each well. Read immediately.

3. PRECAUTIONS

Please read these instructions carefully prior to beginning the assay.

Modifications to the kit components or procedures may result in loss of performance.

4. STORAGE AND STABILITY

Store kit at 4°C immediately upon receipt, apart from the SP Conjugate & Biotinylated Antibody, which should be stored at -20°C.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in sections 9 & 10.

5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
Plasminogen Microplate (12 x 8 well strip)	96 wells	4°C
Plasminogen Standard	1 vial	4°C
10X Diluent M Concentrate	30 mL	4°C
Biotinylated Human Plasminogen Antibody	1 vial	-20°C
100X Streptavidin-Peroxidase Conjugate (SP Conjugate)	80 µL	-20°C
Chromogen Substrate	7 mL	4°C
Stop Solution	11 mL	4°C
20X Wash Buffer Concentrate	2 x 30 mL	4°C
Sealing Tapes	3	N/A

6. MATERIALS REQUIRED, NOT SUPPLIED

These materials are not included in the kit, but will be required to successfully utilize this assay:

- 1 Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 1 μ L to 1 mL volumes.
- Adjustable 1-25 mL pipettes for reagent preparation.
- 100 mL and 1 liter graduated cylinders.
- Absorbent paper.
- Distilled or deionized water.
- Log-log graph paper or computer and software for ELISA data analysis.
- 8 tubes to prepare standard or sample dilutions.

7. LIMITATIONS

- Do not mix or substitute reagents or materials from other kit lots or vendors.

8. TECHNICAL HINTS

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps.
- **This kit is sold based on number of tests. A ‘test’ simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.**

9. REAGENT PREPARATION

Equilibrate all reagents to room temperature (18-25°C) prior to use. Prepare fresh reagents immediately prior to use. If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved.

9.1 1X Diluent M

Dilute the 10X Diluent M Concentrate 1:10 with reagent grade water. Mix gently and thoroughly.

9.2 1X Wash Buffer

Dilute the 20X Wash Buffer Concentrate 1:20 with reagent grade water. Mix gently and thoroughly.

9.3 1X Biotinylated Plasminogen Detector Antibody

9.3.1 The stock Biotinylated Plasminogen Antibody must be diluted with 1X Diluent M according to the label concentration to prepare 1X Biotinylated Plasminogen Antibody for use in the assay procedure. Observe the label for the “X” concentration on the vial of Biotinylated Plasminogen Antibody.

9.3.2 Calculate the necessary amount of 1X Diluent M to dilute the Biotinylated Plasminogen Antibody to prepare a 1X Biotinylated Plasminogen Antibody solution for use in the assay procedure according to how many wells you wish to use and the following calculation:

Number of Wells Strips	Number of Wells	(V _T) Total Volume of 1X Biotinylated Antibody (μL)
4	32	1,760
6	48	2,640
8	64	3,520
10	80	4,400
12	96	5,280

Any remaining solution should be frozen at -20°C.

Where:

C_S = Starting concentration (X) of stock Biotinylated Plasminogen Antibody (variable)

C_F = Final concentration (always = 1X) of 1X Biotinylated Plasminogen Antibody solution for the assay procedure

V_T = Total required volume of 1X Biotinylated Plasminogen Antibody solution for the assay procedure

V_A = Total volume of (X) stock Biotinylated Plasminogen Antibody

V_D = Total volume of 1X Diluent M required to dilute (X) stock Biotinylated Plasminogen Antibody to prepare 1X Biotinylated Plasminogen solution for assay procedures

Calculate the volume of (X) stock Biotinylated Antibody required for the given number of desired wells:

$$(C_F / C_S) \times V_T = V_A$$

Calculate the final volume of 1X Diluent M required to prepare the 1X Biotinylated Plasminogen Antibody:

$$V_T - V_A = V_D$$

Example:

NOTE: This example is for demonstration purposes only. Please remember to check your antibody vial for the actual concentration of antibody provided.

C_S = 50X Biotinylated Plasminogen Antibody stock

C_F = 1X Biotinylated Plasminogen Antibody solution for use in the assay procedure

V_T = 3,520 μ L (8 well strips or 64 wells)

$$(1X/50X) \times 3,520 \mu\text{L} = 70.4 \mu\text{L}$$

$$3,520 \mu\text{L} - 70.4 \mu\text{L} = 3,449.6 \mu\text{L}$$

V_A = 70.4 μ L total volume of (X) stock Biotinylated Plasminogen Antibody required

V_D = 3,449.6 μ L total volume of 1X Diluent M required to dilute the 50X stock Biotinylated Antibody to prepare 1X Biotinylated Plasminogen Antibody solution for assay procedures

9.3.3 First spin the Biotinylated Plasminogen Antibody vial to collect the contents at the bottom.

9.3.4 Add calculated amount V_A of stock Biotinylated Plasminogen Antibody to the calculated amount V_D of 1X Diluent M. Mix gently and thoroughly.

9.4 **1X SP Conjugate**

Spin down the 100X Streptavidin-Peroxidase Conjugate (SP Conjugate) briefly and dilute the desired amount of the conjugate 1:100 with 1X Diluent M.

Any remaining solution should be frozen at -20°C.

10. STANDARD PREPARATIONS

- Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.
- Any remaining standard should be stored at -20°C after reconstitution and used within 30 days.
- This procedure prepares sufficient standard dilutions for duplicate wells.

10.1 Reconstitution of the Plasminogen Standard vial to prepare the 40 ng/mL Plasminogen **Stock Standard**:

10.1.1 First consult the Plasminogen Standard vial to determine the mass of protein in the vial.

10.1.2 Calculate the appropriate volume of 1X Diluent M to add when resuspending the Plasminogen Standard vial to produce a 40 ng/mL Plasminogen **Stock Standard** by using the following equation:

C_S = Starting mass of Plasminogen Standard (see vial label)
(ng)

C_F = 40 ng/mL Plasminogen **Stock Standard** final required concentration

V_D = Required volume of 1X Diluent M for reconstitution (μL)

Calculate total required volume 1X Diluent M for resuspension:
$$(C_S / C_F) \times 1,000 = V_D$$

Example:

NOTE: This example is for demonstration purposes only. Please remember to check your standard vial for the actual amount of standard provided.

C_S = 320 ng of Plasminogen Standard in vial

C_F = 40 ng/mL Plasminogen **Stock Standard** final concentration

V_D = Required volume of 1X Diluent M for reconstitution
$$(320 \text{ ng} / 40 \text{ ng/mL}) \times 1,000 = 8,000 \text{ } \mu\text{L}$$

- 10.1.3 First briefly spin the Plasminogen Standard Vial to collect the contents on the bottom of the tube.
- 10.1.4 Reconstitute the Plasminogen Standard vial by adding the appropriate calculated amount V_D of 1X Diluent M to the vial to generate the 40 ng/mL Plasminogen **Stock Standard**. Mix gently and thoroughly.
- 10.2 Allow the reconstituted 40 ng/mL Plasminogen **Stock Standard** to sit for 10 minutes with gentle agitation prior to making subsequent dilutions.
- 10.3 Add 240 μ L of the **Stock Standard** into tube# **Standard #1**
- 10.4 Label seven tubes #2 – 8
- 10.5 Add 120 μ L of 1X Diluent M to tube #2 – 8.
- 10.6 To prepare **Standard #2**, add 120 μ L of the **Standard #1** into tube #2 and mix gently.
- 10.7 To prepare **Standard #3**, add 120 μ L of the **Standard #2** into tube #3 and mix gently.
- 10.8 Using the table below as a guide, prepare subsequent serial dilutions.
- 10.9 1X Diluent M serves as the zero standard, 0 ng/mL (tube #8).

ASSAY PREPARATION

Standard Dilution Preparation Table

Standard #	Volume to Dilute (μL)	Volume Diluent M (μL)	Total Volume (μL)	Starting Conc. (ng/mL)	Final Conc. (ng/mL)
1	Step 10.1				40.00
2	120	120	240	80.00	20.00
3	120	120	240	40.00	10.00
4	120	120	240	20.00	5.00
5	120	120	240	10.00	2.500
6	120	120	240	5.000	1.250
7	120	120	240	2.500	0.625
8	-	120	120	-	0



11. SAMPLE PREPARATION

11.1 Plasma

Collect plasma using one-tenth volume of 0.1 M sodium citrate as an anticoagulant. (EDTA or Heparin can also be used as an anticoagulant). Centrifuge samples at 3,000 x *g* for 10 minutes. Dilute samples 1:20,000 with 1X Diluent M and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.2 Serum

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 3,000 x *g* for 10 minutes and remove the serum. Dilute samples 1:20,000 with 1X Diluent M and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.3 Saliva

Collect saliva using sample tube. Centrifuge samples at 800 x *g* for 10 minutes and assay. The sample is suggested for use at 1x or within the range of 2x – 10x into Diluent M; however, user should determine optimal dilution factor depending on application needs. Store undiluted samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.4 Urine

Collect urine using sample pot. Centrifuge samples at 800 x *g* for 10 minutes and assay. The sample is suggested for use at 1x or within the range of 2x – 10x into Diluent M; however, user should determine optimal dilution factor depending on application needs. Store undiluted samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.5 **Milk**

Collect milk using sample tube. Centrifuge samples at 800 x *g* for 10 minutes. Milk dilution is suggested at 1:40 in 1X Diluent M then assay or within the range of 5x – 500x; however, the user should determine the optimal dilution factor. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.6 **Cerebrospinal Fluid**

Collect cerebrospinal fluid using sample pot. Centrifuge samples at 3,000 x *g* for 10 minutes. A 1:40 sample dilution is suggested into Diluent M or within the range of 5x – 500x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -80°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

12. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused well plate strips should be returned to the plate packet and stored at 4°C.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

13. ASSAY PROCEDURE

- **Equilibrate all materials and prepared reagents to room temperature (18 - 25°C) prior to use.**
 - **It is recommended to assay all standards, controls and samples in duplicate.**
- 13.1 Prepare all reagents, working standards and samples as instructed. Equilibrate reagents to room temperature before use. The assay is performed at room temperature (18-25°C).
 - 13.2 Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccant inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
 - 13.3 Add 50 µL of Plasminogen Standard or sample per well. Cover wells with a sealing tape and incubate for one hour. Start the timer after the last sample addition.
 - 13.4 Wash five times with 200 µL of 1X Wash Buffer manually. Invert the plate each time and decant the contents; tap it 4-5 times on absorbent paper towel to completely remove the liquid. If using a machine wash six times with 300 µL of 1X Wash Buffer and then invert the plate, decant the contents; tap it 4-5 times on absorbent paper towel to completely remove the liquid.
 - 13.5 Add 50 µL of 1X Biotinylated Plasminogen Antibody to each well and incubate for one hour.
 - 13.6 Wash microplate as described above.
 - 13.7 Add 50 µL of 1X SP Conjugate to each well and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
 - 13.8 Wash microplate as described above.
 - 13.9 Add 50 µL of Chromogen Substrate per well. Gently tap plate to thoroughly coat wells. Break any bubbles that may have

formed. Incubate in ambient light for 30 minutes or until the optimal blue color density develops.

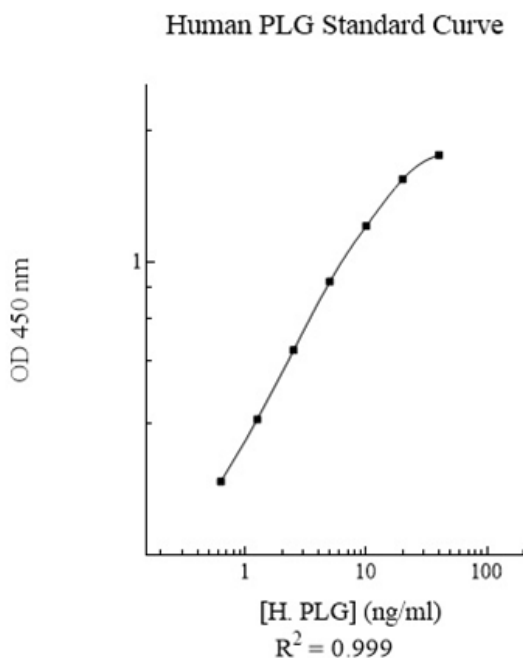
- 13.10 Add 50 μ L of Stop Solution to each well. The color will change from blue to yellow.
- 13.11 Read the absorbance on a microplate reader at a wavelength of 450 nm immediately. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. Please note that some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

14. CALCULATIONS

Calculate the mean value of the triplicate readings for each standard and sample. To generate a Standard Curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance on the y-axis. The best-fit line can be determined by regression analysis using log-log or four-parameter logistic curve-fit. Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.

15. TYPICAL DATA

TYPICAL STANDARD CURVE – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



16. TYPICAL SAMPLE VALUES

SENSITIVITY –

The minimum detectable dose of Plasminogen is typically ~0.23 ng/mL.

SPIKING RECOVERY –

Recovery was determined by spiking one plasma and one serum sample with different human PLG concentrations

Sample	Unspiked Sample (ng/ml)	Spiking Value (ng/ml)	Expected	Observed	Recovery (%)
Plasma	7.103	4.897	12.000	12.183	102
		2.559	9.662	8.832	91
		1.188	8.291	8.252	100
Serum	7.993	4.897	12.890	12.428	96
		2.559	10.552	9.992	95
		1.188	9.181	9.491	103
Average Recovery (%)					98

REFERENCE VALUES

The normal Human plasma and serum levels of Plasminogen is between 70 to 215 µg/mL.

LINEARITY OF DILUTION –

Plasma Dilution	Average % Expected Value
1:10,000	95
1:20,000	99
1:40,000	104

Serum Dilution	Average % Expected Value
1:10,000	96
1:20,000	101
1:40,000	104

PRECISION

	Intra-Assay	Inter-Assay
Average CV (%)	4.3	10.2

17. ASSAY SPECIFICITY

Species	% Cross Reactivity
Canine	None
Bovine	None
Equine	None
Monkey	3
Mouse	None
Rat	None
Swine	None
Rabbit	None
Human	100

Proteins	Cross Reactivity (%)
Human Lp(a)	None
Plasmin	<30%

18. TROUBLESHOOTING

Problem	Cause	Solution
Poor standard curve	Improper standard dilution	Confirm dilutions made correctly
	Standard improperly reconstituted (if applicable)	Briefly spin vial before opening; thoroughly resuspend powder (if applicable)
	Standard degraded	Store sample as recommended
	Curve doesn't fit scale	Try plotting using different scale
Low signal	Incubation time too short	Try overnight incubation at 4°C
	Target present below detection limits of assay	Decrease dilution factor; concentrate samples

RESOURCES

Problem	Cause	Solution
	Precipitate can form in wells upon substrate addition when concentration of target is too high	Increase dilution factor of sample
	Using incompatible sample type (e.g. serum vs. cell extract)	Detection may be reduced or absent in untested sample types
	Sample prepared incorrectly	Ensure proper sample preparation/dilution
Large CV	Bubbles in wells	Ensure no bubbles present prior to reading plate
	All wells not washed equally/thoroughly	Check that all ports of plate washer are unobstructed wash wells as recommended
	Incomplete reagent mixing	Ensure all reagents/master mixes are mixed thoroughly
	Inconsistent pipetting	Use calibrated pipettes and ensure accurate pipetting
	Inconsistent sample preparation or storage	Ensure consistent sample preparation and optimal sample storage conditions (eg. minimize freeze/thaws cycles)

Problem	Cause	Solution
High background/ Low sensitivity	Wells are insufficiently washed	Wash wells as per protocol recommendations
	Contaminated wash buffer	Make fresh wash buffer
	Waiting too long to read plate after adding STOP solution	Read plate immediately after adding STOP solution

RESOURCES

	Improper storage of ELISA kit	Store all reagents as recommended. Please note all reagents may not have identical storage requirements.
	Using incompatible sample type (e.g. Serum vs. cell extract)	Detection may be reduced or absent in untested sample types

19. NOTES

Technical Support

Copyright © 2026 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)