

Version 2c Last updated 3 June 2026

ab325257

Human Fibrinogen

ELISA Kit

For the quantitative measurement of human Fibrinogen in plasma samples.

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

Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	6
9. Reagent Preparation	7
10. Human Fibrinogen Standard Preparation	8
11. Sample Preparation	11
12. Plate Preparation	12
13. Assay Procedure	13
14. Typical Data	15
15. Typical Sample Values	16
16. Assay Specificity	17
17. Species Reactivity	18
18. Troubleshooting	19
19. Notes	21

1. Overview

Human Fibrinogen *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Fibrinogen levels in plasma samples.

A Fibrinogen specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Fibrinogen is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 inversely proportional to the amount of Fibrinogen captured in plate.

Fibrinogen (FBG) is a homodimer of molecular mass 340 kDa, made up of two sets of α , β , γ polypeptide chains, and synthesized in the parenchymal cell of the hepatocyte and in the megakaryocyte. Fibrinogen plays a major role in coagulation, and both elevated and decreased levels have clinical significance. Upon cleavage by thrombin in the initial stages of coagulation activation, Fibrinogen self-assembles to yield a fibrin clot matrix that subsequently is crosslinked by factor XIIIa to form an insoluble network. Fibrinogen also binds to the platelet glycoprotein IIb/IIIa receptor to form bridges between platelets, thus facilitating aggregation. Elevated plasma Fibrinogen has been identified as an independent risk factor for coronary atherosclerosis and ischemic heart disease. Individuals with congenital absence of Fibrinogen, termed afibrinogenemia, have prolonged bleeding times.

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add standard, sample and prepared biotin protein to appropriate wells. Incubate at room temperature.



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



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.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- If applicable, please refer to the current Safety Data Sheet (SDS) provided with this product for safety, handling, and disposal information. The most up to date and current versions are available on our website <https://www.abcam.com/en-us>.

4. Storage and Stability

Store kit at +4°C immediately upon receipt, apart from the SP Conjugate, which should be stored at -20°C. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

5. Limitations

- Assay kit is intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

Item	Quantity	Storage Condition
Human Fibrinogen Microplate	96 wells	4°C
Sealing Tapes	3	N/A
Human Fibrinogen Standard	1 vial	4°C
Biotinylated Human Fibrinogen Protein (2x)	1 vial	4°C
Diluent N Concentrate (10x)	30 mL	4°C
20X Wash Buffer Concentrate	30 mL	4°C
SP Conjugate (100x)	80 µL	-20°C
Chromogen Substrate (1x)	7 mL	4°C
Stop Solution (1x)	11 mL	4°C

7. Materials Required, Not Supplied

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

- Microplate reader is 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.
- 6 tubes to prepare standard or sample dilutions.

8. Technical Hints

- 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.
- Selected components in this kit are supplied in surplus amounts to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Make sure all buffers and solutions are at room temperature before starting the experiment.
- 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.
- Make sure you have the right type of plate for your detection method of choice.
- Make sure the heat block/water bath and microplate reader are switched on before starting the experiment.

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 Diluent N Concentrate (10x)

Dilute the Diluent N Concentrate 10-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved. Store for up to 30 days at 2 - 8°C.

9.2 Biotinylated Human Fibrinogen Protein (2x)

- 9.2.1 Reconstitute the Biotinylated Human Fibrinogen Protein with 3 ml of Diluent N to generate a stock solution. Allow the vial to sit for 10 minutes with gentle agitation prior to dilution.
- 9.2.2 From the stock solution, dilute 2-fold with Diluent N to produce a 1x working solution.
- 9.2.3 Aliquot remaining stock solution to limit repeated freeze-thaw cycles. This solution should be stored at -20°C and used within 5 days.

9.3 20X Wash Buffer Concentrate

Dilute the Wash Buffer Concentrate 20-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved.

9.4 SP Conjugate (100x)

Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 100-fold with Diluent N to produce a 1x solution. The undiluted conjugate should be stored at -20°C.

10. Human Fibrinogen Standard Preparation

- Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.
- Any remaining standard should be aliquoted and stored at -20°C after reconstitution and used within 30 days. Avoid freeze thaw cycles. When needed, the frozen stock solution should be thawed in a water bath at 37°C for 30 minutes or more.
- This procedure prepares sufficient standard dilutions for duplicate wells.

10.1 Reconstitution of the Human Fibrinogen Standard vial to prepare an 80 µg/ml standard stock solution:

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

10.1.2 Calculate the appropriate volume of 1X Diluent N to add when resuspending the Human Fibrinogen Standard vial to produce an 80 µg/ml Human Fibrinogen **Stock Standard Solution** by using the following equation:

C_S = Starting mass of Human Fibrinogen Standard (see vial label)
(µg)

C_F = 80 µg/mL Human Fibrinogen **Stock Standard solution** final
required concentration

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

Calculate total required volume 1X Diluent N 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.

$C_S = 40 \mu\text{g}$ of Human Fibrinogen Standard in vial

$C_F = 80 \mu\text{g/mL}$ Human Fibrinogen **Stock Standard** final concentration

$V_D =$ Required volume of 1X Diluent N for reconstitution

$$(40 \mu\text{g} / 80 \mu\text{g} / \text{mL}) \times 1,000 = 500 \mu\text{L}$$

- 10.1.3 First briefly spin the Human Fibrinogen Standard Vial to collect the contents on the bottom of the tube.
- 10.1.4 Reconstitute the Human Fibrinogen Standard vial by adding the appropriate calculated amount V_D of 1X Diluent N to the vial to generate the $80 \mu\text{g/mL}$ Human Fibrinogen **Stock Standard solution**. Mix gently and thoroughly.
- 10.2 Allow the reconstituted $80 \mu\text{g/mL}$ Human Fibrinogen **Stock Standard solution** to sit for 10 minutes with gentle agitation prior to making subsequent dilutions
- 10.3 Label eight tubes #2-6.
- 10.4 Prepare the $40 \mu\text{g/mL}$ **Standard #2** by adding $120 \mu\text{L}$ of the reconstituted $80 \mu\text{g/mL}$ Human Fibrinogen **Stock Standard solution** to $120 \mu\text{L}$ of 1X Diluent N and mix thoroughly and gently.
- 10.5 Add $120 \mu\text{L}$ of 1X Diluent N to tubes #3 – 6.
- 10.6 To prepare **Standard #3**, add $120 \mu\text{L}$ of the **Standard #2** into tube #3 and mix gently.
- 10.7 To prepare **Standard #4**, add $120 \mu\text{L}$ of the **Standard #3** into tube #4 and mix gently.
- 10.8 Using the table below as a guide, prepare subsequent serial dilutions.
- 10.9 1X Diluent N serves as the zero standard, $0 \mu\text{g/mL}$ (tube #6).

Standard Dilution Preparation Table

Standard #	Volume to Dilute (μL)	Volume Diluent N (μL)	Total Volume (μL)	Starting Conc. (μg/mL)	Final Conc. (μg/mL)
1	Step 10.1				80
2	120	120	240	80	40
3	120	120	240	40	20
4	120	120	240	20	10
5	120	120	240	10	5
6	-	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 3000 x g for 10 minutes and collect plasma. A 250-fold sample dilution is suggested into Diluent N; however, users should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

ΔNote: Applicable samples may also include biofluids, cell culture, and tissue homogenates. If necessary, users should determine optimal dilution factor depending on application needs.

Refer to Dilution Guidelines for further instruction.

Guidelines for Dilutions of 100-fold or Greater <i>(for reference only; please follow the insert for specific dilution suggested)</i>	
100x	10000x
4 µl sample + 396 µl buffer (100X) = 100-fold dilution <i>Assuming the needed volume is less than or equal to 400 µl</i>	A) 4 µl sample + 396 µl buffer (100X) B) 4 µl of A + 396 µl buffer (100X) = 10000-fold dilution <i>Assuming the needed volume is less than or equal to 400 µl</i>
1000x	100000x
A) 4 µl sample + 396 µl buffer (100X) B) 24 µl of A + 216 µl buffer (10X) = 1000-fold dilution <i>Assuming the needed volume is less than or equal to 240 µl</i>	A) 4 µl sample + 396 µl buffer (100X) B) 4 µl of A + 396 µl buffer (100X) C) 24 µl of A + 216 µl buffer (10X) = 100000-fold dilution <i>Assuming the needed volume is less than or equal to 240 µl</i>

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.

13. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
 - We recommend that you assay all standards, controls and samples in duplicate.
- 13.1** Prepare all reagents, working standards, and samples as directed in the previous sections. The assay is performed at room temperature (20-25°C).
 - 13.2** Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
 - 13.3** Add 25 µL of Fibrinogen Standard or sample to each well and immediately add 25 µL of Biotinylated Human Fibrinogen Protein to each well (on top of standard or sample). Gently tap plate to ensure thorough mixing. Cover wells with a sealing tape and incubate for two hours. Start the timer after the last sample addition.
 - 13.4** Wash the microplate manually or automatically using a microplate washer. Invert the plate and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If washing manually, wash five times with 200 µl of Wash Buffer per well. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a microplate washer, wash six times with 300 µl of Wash Buffer per well; invert the plate and hit 4-5 times on absorbent material to completely remove the liquid.
 - 13.5** Add 50 µl of SP Conjugate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
 - 13.6** Wash the microplate as described above.
 - 13.7** Add 50 µl of Chromogen Substrate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate in ambient light for 12 minutes or until the optimal blue color density develops.
 - 13.8** Add 50 µl of Stop Solution to each well. The color will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.

13.9 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 low concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

14. Typical Data

Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.

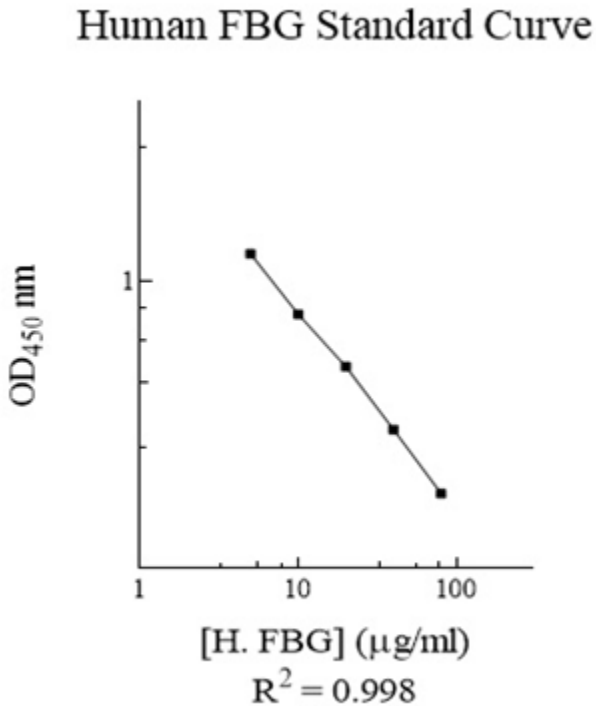


Figure 1. Example of Fibrinogen standard curve. The standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

15. Typical Sample Values

SENSITIVITY –

The minimum detectable dose (MDD) of Fibrinogen as calculated by 2 standard deviations from the mean of a zero standard was established to be 2.9 µg/ml.

PRECISION –

Intra-assay precision was determined by testing replicates of three plasma samples twenty times in one assay.

Inter-assay precision was determined by testing three plasma samples in twenty assays.

	Intra-assay Precision	Inter-Assay Precision
CV (%)	4.7	10.8

SPIKING RECOVERY –

Recovery was determined by spiking two plasma samples with different FBG concentrations.

Sample	Unspiked Sample (µg/mL)	Spiking Value (µg/mL)	Expected	Observed	Recovery (%)
1	6.599	7.871	14.470	13.172	91%
		3.631	10.230	10.629	104%
		1.867	8.466	7.626	90%
2	6.245	7.871	14.116	13.493	96%
		3.631	9.876	8.917	90%
		1.867	8.112	7.533	93%
Average Recovery (%)					94%

LINEARITY OF DILUTION -

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Plasma samples were serially-diluted to test for linearity.

Average Percentage of Expected Value (%)	
Sample Dilution	Plasma
125x	90%
250x	98%
500x	110%

16. Assay Specificity

This kit recognizes Fibrinogen in plasma.

REFERENCE VALUE –

Normal human plasma Fibrinogen concentration has been reported ranging approximately 1.5 to 4 mg/mL. Plasma samples from healthy adults were tested (n=10). On average, human FBG level was 2.6 mg/ml.

17. Species Reactivity

Species	% Cross Reactivity
Canine	None
Bovine	None
Monkey	None
Mouse	None
Rat	None
Swine	None
Rabbit	None

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
	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
	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)