

Mouse Fibrinogen ELISA Kit ab108844

[4 References](#) [2 Images](#)

Overview

Product name Mouse Fibrinogen ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				5.2%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				10%

Sample type Plasma

Assay type Competitive

Sensitivity = 0.6 µg/ml

Range 0.625 µg/ml - 20 µg/ml

Recovery 99 %

Assay time 3h 0m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview

Abcam's Fibrinogen Mouse *in vitro* competitive ELISA (enzyme-linked immunosorbent assay) kit is designed for the quantitative measurement of mouse fibrinogen in plasma.

A Fibrinogen specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a 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.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Platform Microplate

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
2X Biotinylated Mouse Fibrinogen (Lyophilized)	1 vial
20X Wash Buffer Concentrate	1 x 30ml
Chromogen Substrate	1 x 7ml
Fibrinogen Microplate (12 x 8 well strips)	1 unit
Fibrinogen Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function Fibrinogen has a double function: yielding monomers that polymerize into fibrin and acting as a cofactor in platelet aggregation.

Tissue specificity Plasma.

Involvement in disease Defects in FGA are a cause of congenital afibrinogenemia (CAFBN) [MIM:202400]. This is a rare autosomal recessive disorder characterized by bleeding that varies from mild to severe and by complete absence or extremely low levels of plasma and platelet fibrinogen. Note=The majority of cases of afibrinogenemia are due to truncating mutations. Variations in position Arg-35 (the site of cleavage of fibrinopeptide a by thrombin) leads to alpha-dysfibrinogenemias. Defects in FGA are a cause of amyloidosis type 8 (AMYL8) [MIM:105200]; also known as systemic non-neuropathic amyloidosis or Ostertag-type amyloidosis. AMYL8 is a hereditary generalized amyloidosis due to deposition of apolipoprotein A1, fibrinogen and lysozyme amyloids. Viscera are particularly affected. There is no involvement of the nervous system. Clinical features include renal amyloidosis resulting in nephrotic syndrome, arterial hypertension, hepatosplenomegaly, cholestasis, petechial skin rash.

Sequence similarities Contains 1 fibrinogen C-terminal domain.

Domain A long coiled coil structure formed by 3 polypeptide chains connects the central nodule to the C-terminal domains (distal nodules). The long C-terminal ends of the alpha chains fold back, contributing a fourth strand to the coiled coil structure.

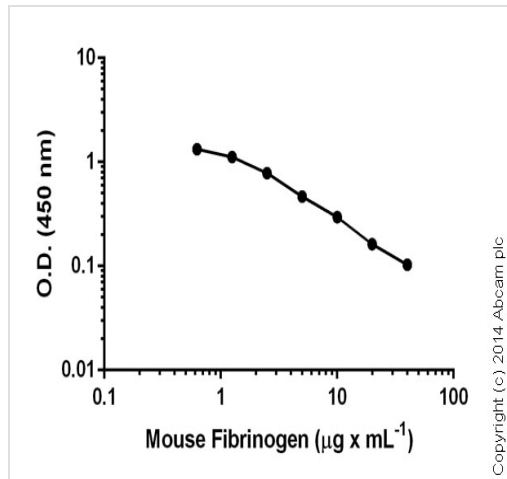
Post-translational modifications The alpha chain is not glycosylated. Forms F13A-mediated cross-links between a glutamine and the epsilon-amino group of a lysine residue, forming fibronectin-fibrinogen heteropolymers.

About one-third of the alpha chains in the molecules in blood were found to be phosphorylated. Conversion of fibrinogen to fibrin is triggered by thrombin, which cleaves fibrinopeptides A and B from alpha and beta chains, and thus exposes the N-terminal polymerization sites responsible for the formation of the soft clot. The soft clot is converted into the hard clot by factor XIIIa which catalyzes the epsilon-(gamma-glutamyl)lysine cross-linking between gamma chains (stronger) and between alpha chains (weaker) of different monomers. Phosphorylation sites are present in the extracellular medium.

Cellular localization

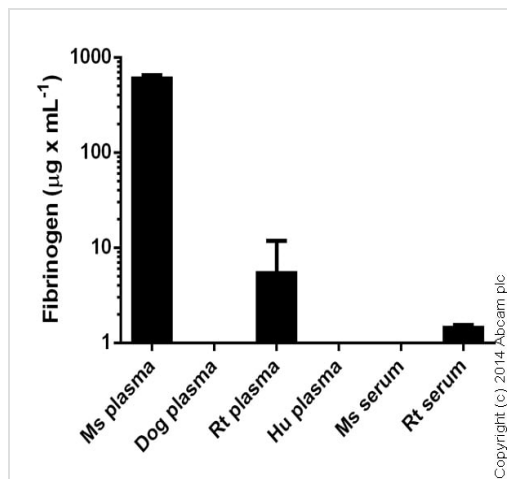
Secreted.

Images



Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Competitive ELISA - Fibrinogen Mouse ELISA Kit (ab108844)



Fibrinogen measured in biological fluids showing quantity (μg) per mL of tested sample. Samples diluted 1-500 fold.

Competitive ELISA - Fibrinogen Mouse ELISA Kit (ab108844)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors