

Product datasheet

Human Fibrinogen ELISA Kit ab108841

★★★★★ 1 Abreviews 7 References 1 Image

Overview

Product name	Human Fibrinogen ELISA Kit										
Detection method	Colorimetric										
Precision	Intra-assay										
	<table border="1"> <thead> <tr> <th>Sample</th> <th>n</th> <th>Mean</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>4%</td> </tr> </tbody> </table>	Sample	n	Mean	SD	CV%	Overall				4%
Sample	n	Mean	SD	CV%							
Overall				4%							
	Inter-assay										
	<table border="1"> <thead> <tr> <th>Sample</th> <th>n</th> <th>Mean</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>8.6%</td> </tr> </tbody> </table>	Sample	n	Mean	SD	CV%	Overall				8.6%
Sample	n	Mean	SD	CV%							
Overall				8.6%							
Sample type	Cell culture supernatant, Saliva, Milk, Urine										
Assay type	Sandwich (quantitative)										
Sensitivity	= 1 ng/ml										
Range	1.25 ng/ml - 80 ng/ml										
Recovery	96 %										
Assay time	4h 00m										
Assay duration	Multiple steps standard assay										
Species reactivity	Reacts with: Human										
Product overview	Human Fibrinogen ELISA kit is designed for the quantitative measurement of fibrinogen concentrations in cell culture media, saliva, milk and urine 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 a Fibrinogen 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 Fibrinogen captured in plate.

Get better reproducibility in only 90 minutes with Human Fibrinogen ELISA Kit ([ab208036](#)) from our SimpleStep ELISA[®] range.

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
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Fibrinogen Antibody	1 x 120µl
Chromogen Substrate	1 x 8ml
Fibrinogen Microplate (12 x 8 well strips)	1 unit
Fibrinogen Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 12ml

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.

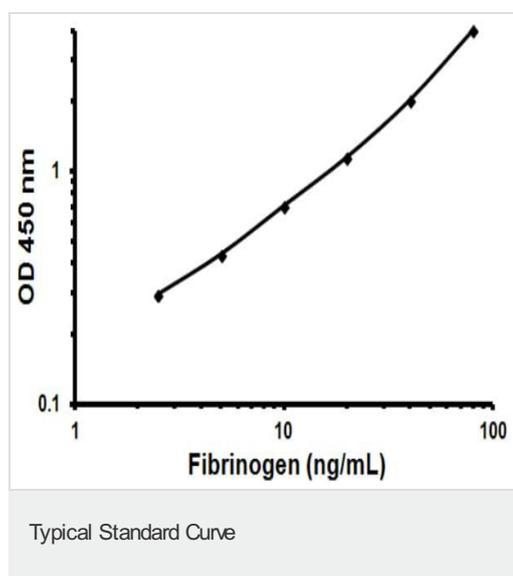
Conversion of fibrinogen to fibrin is triggered by thrombin, which cleaves fibrinopeptides A and B from alpha and beta chains, and thus exposes the N-terminal polymerization sites responsible for the formation of the soft clot. The soft clot is converted into the hard clot by factor XIIIa which catalyzes 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



Representative Standard Curve using ab108841.

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