

Human Tissue Factor Pathway Inhibitor ELISA Kit (TFPI) ab108904

6 References 2 Images

Overview

Product name	Human Tissue Factor Pathway Inhibitor ELISA Kit (TFPI)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			5%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			9.9%
Sample type	Cell culture supernatant, Milk, Serum, Plasma, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	> 0.081 ng/ml			
Range	0.625 ng/ml - 5 ng/ml			
Recovery	98 %			
Assay time	5h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Tissue Factor Pathway Inhibitor (TFPI) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Tissue Factor Pathway Inhibitor concentrations in plasma, serum, milk, cerebrospinal fluid and cell culture samples.			

A Tissue Factor Pathway Inhibitor specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Tissue Factor Pathway Inhibitor 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 Tissue Factor Pathway Inhibitor captured in plate.

Get results in 90 minutes with Human Tissue Factor Pathway Inhibitor ELISA Kit ([ab188392](#)) 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 Tissue Factor Pathway Inhibitor Antibody	1 x 120µl
Chromogen Substrate	1 x 8ml
Sealing Tapes	3 units
Stop Solution	1 x 12ml
Tissue Factor Pathway Inhibitor Microplate (12 x 8 well strips)	1 x 96 units
Tissue Factor Pathway Inhibitor Standard	2 vials

Function Inhibits factor X (X(a)) directly and, in a Xa-dependent way, inhibits VIIa/tissue factor activity, presumably by forming a quaternary Xa/LACI/VIIa/TF complex. It possesses an antithrombotic action and also the ability to associate with lipoproteins in plasma.

Tissue specificity Mostly in endothelial cells.

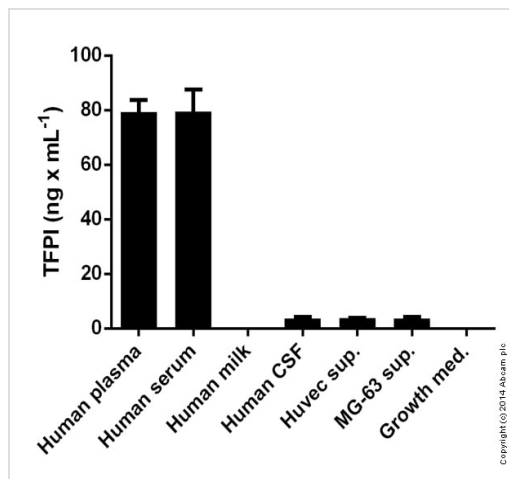
Sequence similarities Contains 3 BPTI/Kunitz inhibitor domains.

Domain This inhibitor contains three inhibitory domains. The first domain interacts with VIIa and TF, the second one with Xa.

Post-translational modifications O-glycosylated.

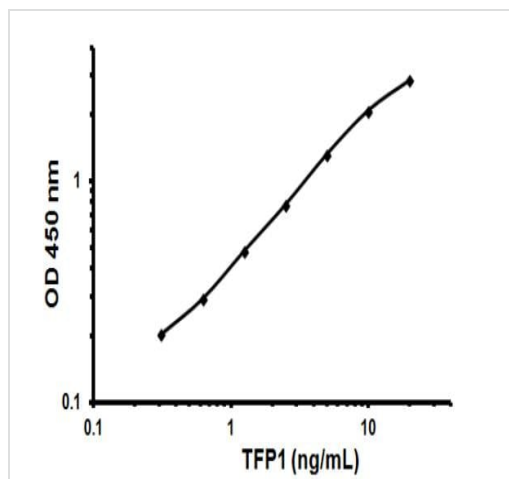
Cellular localization Secreted.

Images



TFPI measured in various samples showing quantity (ng) per mL of tested sample

Sandwich ELISA - ab108904 Tissue Factor Pathway Inhibitor (TFPI) Human ELISA Kit



Representative Standard Curve using ab108904.

Typical Standard Curve

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