

Human Tissue Factor ELISA Kit ab220653

リコンビナント SimpleStep ELISA

4 References [画像数 7](#)

製品の概要

製品名 Human Tissue Factor ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall	8			4.9%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	3			6.9%

サンプルの種類 Cell culture supernatant, Saliva, Urine, Cell culture extracts, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 3.6 pg/ml

検出範囲 15.63 pg/ml - 1000 pg/ml

添加回収試験 特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	110	105% - 116%
Saliva	86	83% - 87%
Urine	83	80% - 85%
Cell culture extracts	105	100% - 107%
Hep Plasma	109	107% - 113%
EDTA Plasma	115	111% - 118%

サンプルの種類	平均 %	測定範囲
Cit plasma	104	99% - 113%

全工程の試験時間

1h 30m

ステップ

One step assay

種交差性

交差種: Human

非交差種: Cow

製品の概要

Human Tissue Factor ELISA Kit (ab220653) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Tissue Factor protein in cell culture extracts, cell culture supernatant, cit plasma, edta plasma, hep plasma, saliva, and urine. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Tissue Factor with 3.6 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

特記事項

Tissue factor (TF), also known as Coagulation factor III, F3 thromboplastin, and CD142 is a 263-amino acid single pass type I membrane protein encoded by the gene F3. TF is a 47 kDa single pass type I membrane protein consisting of an extracellular, transmembrane, and cytoplasmic domain. Tissue factor's main role is in blood coagulation. The protein acts as a receptor for coagulation factor VII, which then forms the active complex FVIIa. This complex proteolytically activates downstream factors, including coagulation Factor IX and X, which leads to the formation of thrombin and fibrin clot formation. Tissue Factor is in a class of proteins known as Cytokine Receptor class II family. This family is activated by cytokines, and play a role in angiogenesis and apoptosis.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

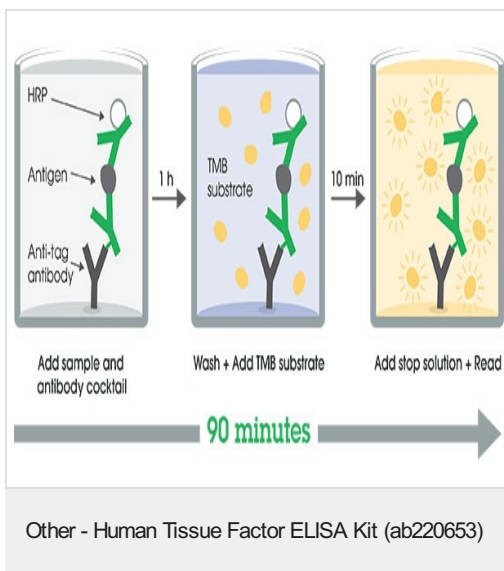
内容	1 x 96 tests
10X Human Tissue Factor Capture Antibody	1 x 600ml
10X Human Tissue Factor Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BI	1 x 6ml
Human Tissue Factor Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

機能 Initiates blood coagulation by forming a complex with circulating factor VII or VIIa. The [TF:VIIa] complex activates factors IX or X by specific limited proteolysis. TF plays a role in normal hemostasis by initiating the cell-surface assembly and propagation of the coagulation protease cascade.

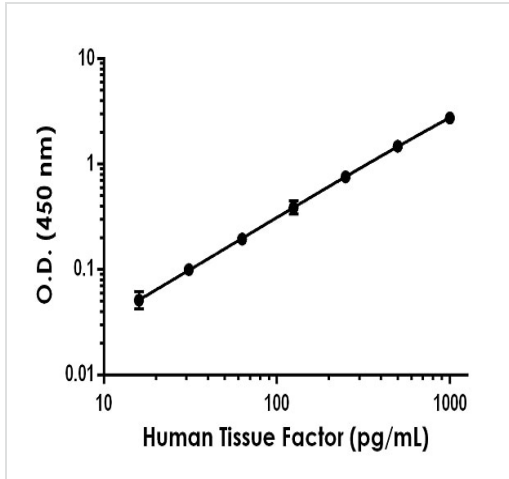
配列類似性 Belongs to the tissue factor family.

細胞内局在 Membrane.

画像

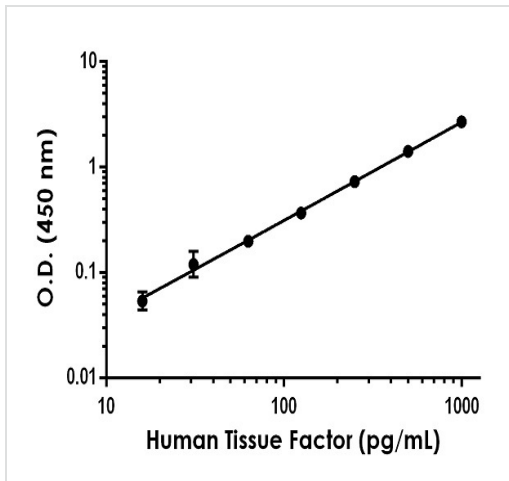


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



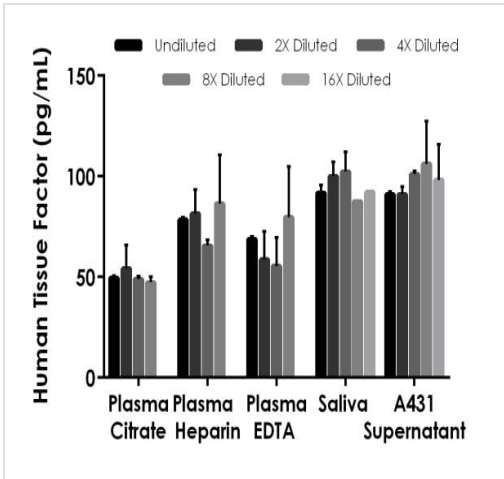
Background-subtracted data values (mean +/- SD) are graphed.

Example of human Tissue Factor standard curve in Sample Diluent NS.



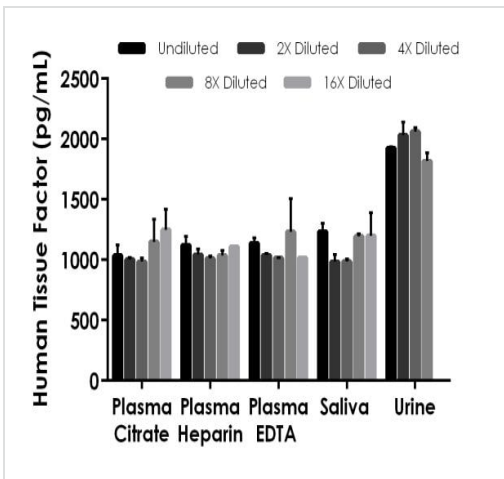
Background-subtracted data values (mean +/- SD) are graphed.

Example of human Tissue Factor standard curve in 1X Cell Extraction Buffer PTR.



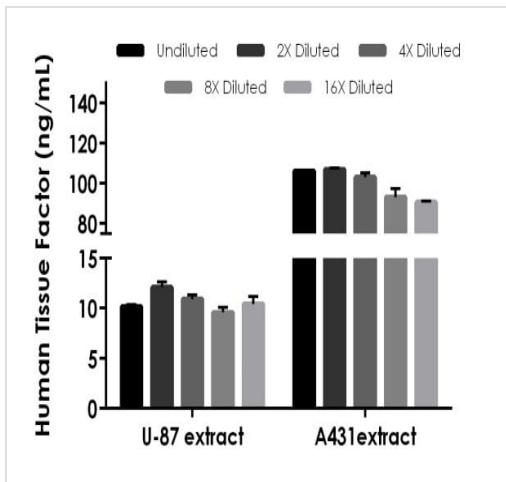
Interpolated concentrations of native Tissue Factor in human plasma and A431 cell culture supernatant samples.

The concentrations of Tissue factor were measured in duplicates, interpolated from the Tissue Factor standard curves and corrected for sample dilution. Undiluted samples are as follows: plasma (citrate) 100%, plasma (heparin) 100%, plasma (EDTA) 100%, saliva 50% and A431 cell culture supernatant 100%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Tissue Factor concentration was determined to be 50 pg/mL in plasma (citrate), 78 pg/mL in plasma (heparin), 65.7 pg/mL in plasma (EDTA), 98 pg/mL in saliva) and 97.5 pg/mL in A431 cell culture supernatant.



Interpolated concentrations of spiked Tissue Factor in human plasma, saliva, and urine samples.

The concentrations of Tissue Factor were measured in duplicates, interpolated from the Tissue Factor standard curves and corrected for sample dilution. Undiluted samples are as follows: 50%, plasma (citrate) 50%, plasma (heparin) 50%, saliva 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Tissue Factor concentration was determined to be 1,000 pg/mL in neat plasma (citrate), plasma (heparin), plasma (EDTA), saliva and 2,000 pg/mL in neat urine.



Interpolated concentrations of native Tissue Factor in U-87 and A431 cellular extract samples based on a 1,000 µg/mL extract load.

The concentrations of Tissue Factor were measured in duplicate and interpolated from the Tissue Factor standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Tissue Factor concentration was determined to be 10.4 ng/mL in U-87 cell extract and 100 ng/mL in A431 cell extract.

Powered by recombinant antibodies

Research with confidence
Consistent and reproducible results

Long-term and scalable supply
Recombinant technology

Success from the first experiment
Confirmed specificity

Ethical standards compliant
Animal-free production

Sandwich ELISA - Human Tissue Factor ELISA Kit (ab220653)

To learn more about the advantages of recombinant antibodies see [here](#).

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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