

Tissue Factor Activity Assay Kit (Human, Colorimetric) ab108906

[16 References](#) [画像数 4](#)

製品の概要

製品名	Tissue Factor Activity Assay Kit (Human, Colorimetric)
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Plasma, Tissue, Cell Lysate
アッセイタイプ	Enzyme activity (quantitative)
全工程の試験時間	1h 00m
種交差性	交差種: Human
製品の概要	Tissue Factor Activity Assay Kit (Human, Colorimetric) ab108906 is used to determine Human Tissue Factor activity in plasma, tissue and cell culture supernatants.

The tissue factor activity assay protocol measures the ability of lipoprotein TF/FVIIa to activate factor X (FX) to factor Xa. The amidolytic activity of the TF/FVIIa complex is quantitated by the amount of FXa produced using a highly specific FXa substrate releasing a yellow para-nitroaniline (pNA) chromophore. The change in absorbance of the pNA at 405 nm is directly proportional to the TF enzymatic activity.

Tissue factor activity assay protocol summary:

- add reaction mix to wells
- add standards and samples
- incubate for 30 min
- add FXa substrate
- analyze with a microplate reader every 5 min for 25 min

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

特記事項

Previously called Tissue Factor Human Chromogenic Activity Assay Kit.

The transmembrane protein Tissue factor (TF) is the physiologic trigger of coagulation in normal hemostasis. Tissue factor binds and allosterically activates factor VII (FVII). The TF-FVIIa complex cleaves factor IX and X, leading to thrombin generation. Tissue factor markedly enhances the ability of FVIIa to cleave both macromolecule and small peptidyl substrates.

試験プラットフォーム

Microplate reader

製品の特性

保存方法

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

内容	100 tests
Assay Diluent	1 x 20ml
FXa Substrate (Lyophilized)	2 vials
Human FVII (Lyophilized)	1 vial
Human FX (Lyophilized)	1 vial
rh Tissue Factor Standard (lipoprotein) (Lyophilized)	1 vial
Sample Diluent	1 x 11ml
Sealing Tapes	3 units
Tissue Factor Microplate (12 strips of 8 wells)	1 unit

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.

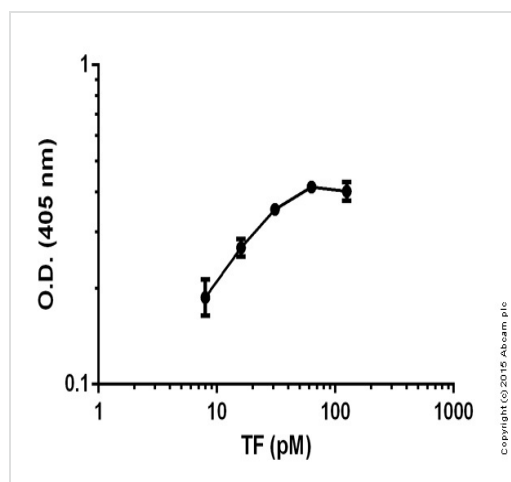
配列類似性

Belongs to the tissue factor family.

細胞内局在

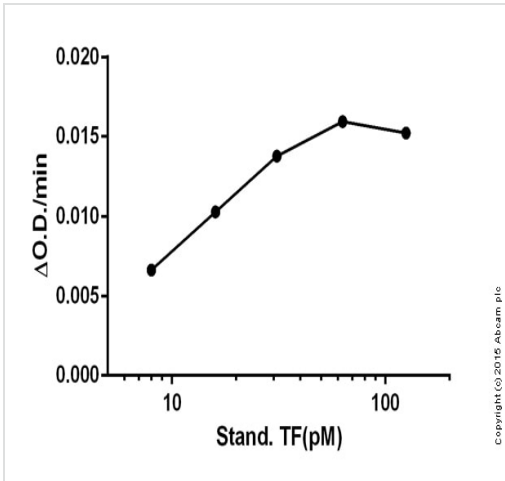
Membrane.

画像



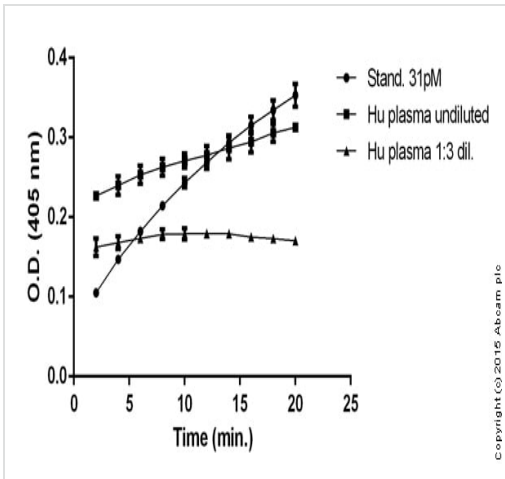
Colorimetric standard curve: mean of duplicates (+/- SD)

Functional studies - Tissue Factor Human
Chromogenic Activity Assay Kit



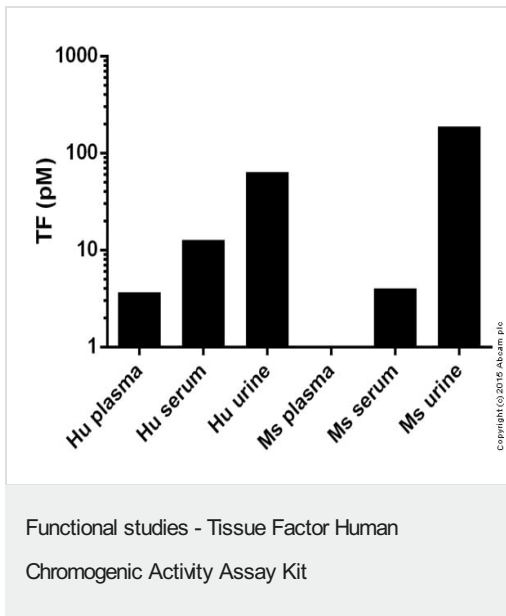
Standard curve showing concentration (pM) on X-axis and increase of O.D. per min on Y-axis (time frame: 20 min.)

Functional studies - Tissue Factor Human
Chromogenic Activity Assay Kit



Time course showing O.D.'s of one standard point and one sample (undiluted vs. diluted)

Functional studies - Tissue Factor Human
Chromogenic Activity Assay Kit



Tissue Factor measured in various biologicals showing average concentrations

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