

Human Tissue Type Plasminogen Activator ELISA Kit (TPA) ab108914

6 References 画像数 3

製品の概要

製品名	Human Tissue Type Plasminogen Activator ELISA Kit (TPA)																				
検出方法	Colorimetric																				
再現性	Intra-Assay (同時再現性) <table><tr><th>サンプル</th><th>N</th><th>平均値</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>2.5%</td></tr></table> Inter-Assay (日差再現性) <table><tr><th>サンプル</th><th>N</th><th>平均値</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td></td><td></td><td></td><td>8.6%</td></tr></table>	サンプル	N	平均値	SD	CV%	Overall				2.5%	サンプル	N	平均値	SD	CV%	Overall				8.6%
サンプル	N	平均値	SD	CV%																	
Overall				2.5%																	
サンプル	N	平均値	SD	CV%																	
Overall				8.6%																	
サンプルの種類	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Tissue Extracts, Cell Lysate, Cerebral Spinal Fluid																				
アッセイタイプ	Sandwich (quantitative)																				
検出感度	= 938 pg/ml																				
検出範囲	60 pg/ml - 500 pg/ml																				
添加回収試験	98 %																				
全工程の試験時間	4h 0m																				
ステップ	Multiple steps standard assay																				
種交差性	交差種: Human																				
製品の概要	Abcam's Tissue Type Plasminogen Activator (TPA) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Tissue Type Plasminogen Activator in plasma, serum, urine, saliva, milk, cell culture supernatants and tissue extracts.																				

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

Get results in 90 minutes with Human Tissue Plasminogen Activator ELISA Kit (**ab190812**) 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.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	1 x 96 tests
100X Biotinylated Human TPA Antibody	1 x 60µl
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml
TPA Microplate (12 x 8 well strips)	1 unit
TPA Standard	1 vial

機能

Converts the abundant, but inactive, zymogen plasminogen to plasmin by hydrolyzing a single Arg-Val bond in plasminogen. By controlling plasmin-mediated proteolysis, it plays an important role in tissue remodeling and degradation, in cell migration and many other physiopathological events. Play a direct role in facilitating neuronal migration.

組織特異性

Synthesized in numerous tissues (including tumors) and secreted into most extracellular body fluids, such as plasma, uterine fluid, saliva, gingival crevicular fluid, tears, seminal fluid, and milk.

関連疾患

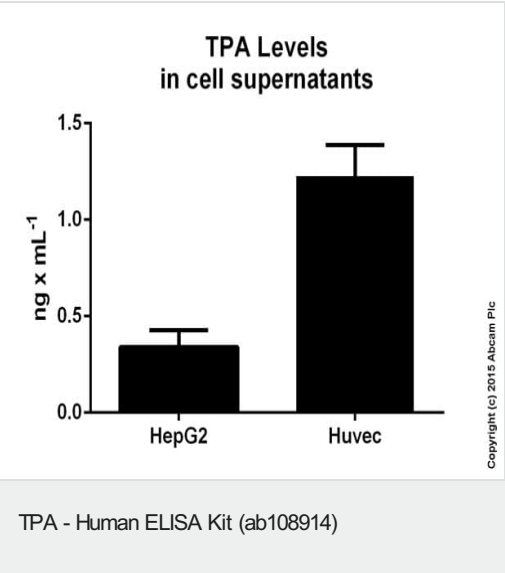
Note=Increased activity of TPA results in increased fibrinolysis of fibrin blood clots that is associated with excessive bleeding. Defective release of TPA results in hypofibrinolysis that can lead to thrombosis or embolism.

配列類似性

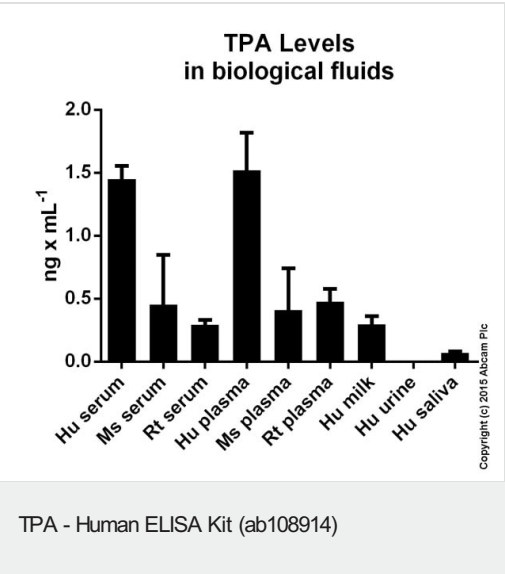
Belongs to the peptidase S1 family.
Contains 1 EGF-like domain.
Contains 1 fibronectin type-I domain.
Contains 2 kringle domains.
Contains 1 peptidase S1 domain.

ドメイン	<p>Both FN1 and one of the kringle domains are required for binding to fibrin.</p> <p>Both FN1 and EGF-like domains are important for binding to LRP1.</p> <p>The FN1 domain mediates binding to annexin A2.</p> <p>The second kringle domain is implicated in binding to cytokeratin-8 and to the endothelial cell surface binding site.</p>
翻訳後修飾	<p>The single chain, almost fully active enzyme, can be further processed into a two-chain fully active form by a cleavage after Arg-310 catalyzed by plasmin, tissue kallikrein or factor Xa.</p> <p>Differential cell-specific N-linked glycosylation gives rise to two glycoforms, type I (glycosylated at Asn-219) and type II (not glycosylated at Asn-219). The single chain type I glycoform is less readily converted into the two-chain form by plasmin, and the two-chain type I glycoform has a lower activity than the two-chain type II glycoform in the presence of fibrin.</p> <p>N-glycosylation of Asn-152; the bound oligomannosidic glycan is involved in the interaction with the mannose receptor.</p> <p>Characterization of O-linked glycan was studied in Bowes melanoma cell line.</p>
細胞内局在	Secreted > extracellular space.

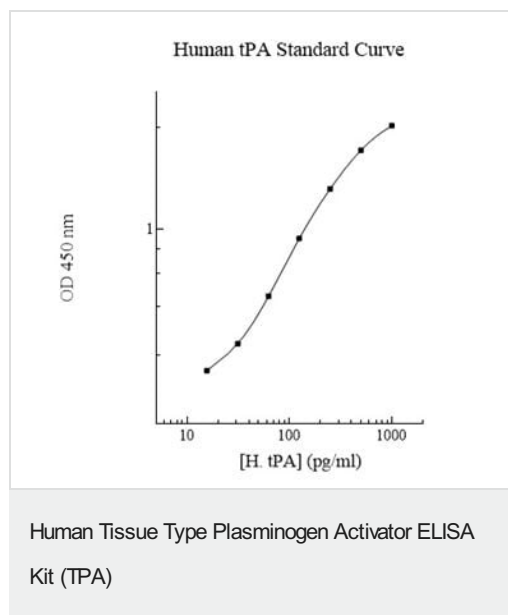
画像



Tissue Type Plasminogen Activator measured in biological fluids showing quantity (ng) per mL of tested sample. Samples diluted 2-10 fold.



Tissue Type Plasminogen Activator measured in biological fluids showing quantity (ng) per mL of tested sample. Human samples diluted 3-27 fold. Rat and mouse samples diluted 1-10 fold.



Representative Standard Curve Using ab108914.

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