

Rat Tissue Plasminogen Activator ELISA Kit (TPA) ab198510

3 References [画像数 2](#)

医薬用外劇物

製品の概要

製品名 Rat Tissue Plasminogen Activator ELISA Kit (TPA)

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%
Known Conc	20	0.8ng/ml	0.034	4.28%
Known Conc	20	4.25ng/ml	0.225	5.29%
Known Conc	20	13.2ng/ml	0.553	4.2%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Known Conc	10	1.37ng/ml	0.06	4.71%
Known Conc	10	4.8ng/ml	0.34	7.13%
Known Conc	10	13.3ng/ml	0.75	5.65%

サンプルの種類 Plasma, Other biological fluids

アッセイタイプ Quantitative

検出感度 0.043 ng/ml

検出範囲 0.05 ng/ml - 50 ng/ml

添加回収試験 99 %

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
Spike	99	93% - 107%

ステップ Multiple steps standard assay

種交差性

交差種: Rat

非交差種: Sheep, Rabbit, Horse, Dog, Human

製品の概要

Abcam's Tissue Plasminogen Activator (TPA) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of active tissue Plasminogen Activator (TPA) in rat plasma and other biological fluids.

Functionally active rat TPA will form a covalent complex with the biotinylated Human PAI-1, which is bound to the avidin on the plate. Only free active TPA will react with the PAI-1 bound to the plate. After appropriate washing steps, polyclonal anti-TPA primary antibody binds to the captured TPA. Excess antibody is washed away and bound primary antibody is then bound by the secondary antibody conjugated to horseradish peroxidase. Following an additional washing step, TMB substrate is used for color development at 450nm. The amount of color development is directly proportional to the concentration of active TPA in the sample.

試験プラットフォーム

Microplate

製品の特性**保存方法**

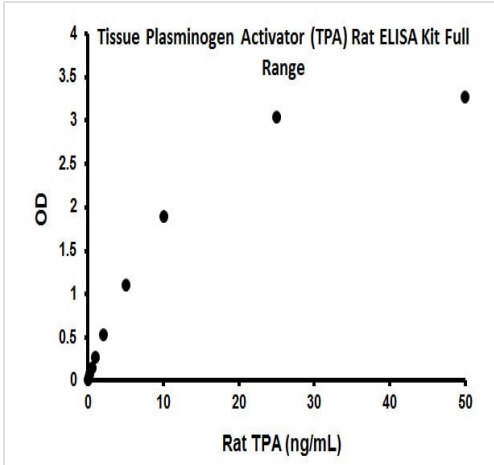
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X TBS buffer pH 7.4	1 x 5ml
10X Wash Buffer	1 x 50ml
HRP secondary antibody	1 x 10µl
Anti-TPA primary antibody	1 vial
Avidin pre-coated 96 well microplate (12 x 8 well strips)	1 x 96 tests
Biotinylated Human PAI-1	1 vial
Rat TPA activity standard	1 vial
TMB Substrate Solution	1 x 10ml

細胞内局在

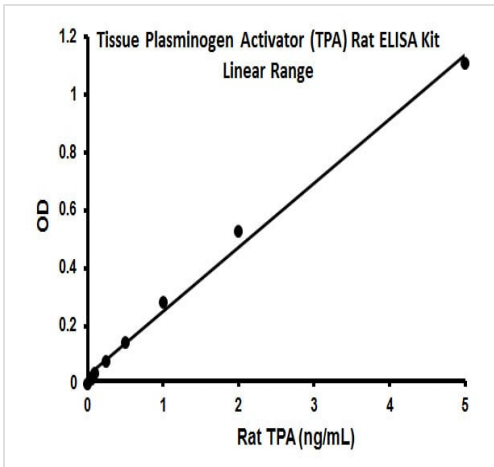
Secreted

画像



Full range standard curve using ab198510 Tissue Plasminogen Activator (TPA) Rat ELISA Kit.

Standard Curve



Linear range standard curve using ab198510 Tissue Plasminogen Activator (TPA) Rat ELISA Kit.

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.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