

Mouse uPA Total ELISA Kit (URK) ab198512

3 References 画像数 2

医薬用外劇物

製品の概要

製品名 Mouse uPA Total ELISA Kit (URK)

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 | 0.06ng/ml | 0.005 | 8.12% |
| Known Conc | 20 | 0.3ng/ml | 0.019 | 6.18% |
| Known Conc | 20 | 3.01ng/ml | 0.112 | 3.73% |

Inter-Assay (日差再現性)

| サンプル | N | 平均値 | SD | CV% |
|------------|----|-----------|-------|-------|
| Known Conc | 10 | 0.09ng/ml | 0.008 | 8.52% |
| Known Conc | 10 | 0.25ng/ml | 0.019 | 7.47% |
| Known Conc | 10 | 2.66ng/ml | 0.154 | 5.79% |

サンプルの種類 Plasma

アッセイタイプ Quantitative

検出感度 0.00998 ng/ml

検出範囲 0.01 ng/ml - 10 ng/ml

添加回収試験 97.75 %

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 | = 97.7 | 89.3% - 106% |

ステップ Multiple steps standard assay

種交差性

交差種: Mouse

非交差種: Sheep

製品の概要

Abcam's uPA Total (URK) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of total urokinase plasminogen activator antigen in mouse plasma.

Mouse uPA will bind to the capture antibody coated on the microtiter plate. Free and complexed enzyme will react with the capture antibody on the plate. After appropriate washing steps, polyclonal anti-murine uPA primary antibody binds to the captured enzyme. Excess antibody is washed away and bound polyclonal antibody is then bound by the secondary antibody conjugated to horseradish peroxidase. TMB substrate is used for color development at 450nm. A standard calibration curve is prepared along with the samples to be measured using dilutions of uPA.

試験プラットフォーム

Microplate

製品の特性

保存方法

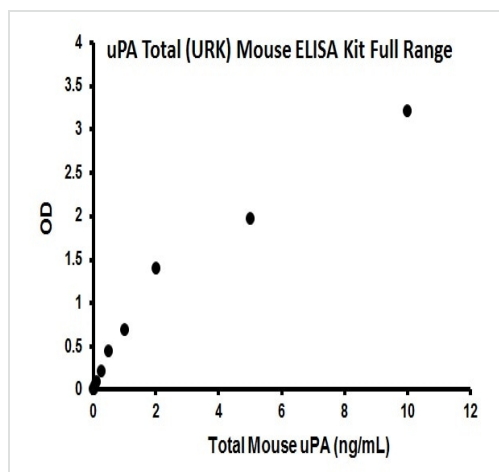
Store at +4°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|---|--------------|
| 10X Wash Buffer | 1 x 50ml |
| Anti-mouse uPA primary antibody | 1 vial |
| Anti-rabbit horseradish peroxidase secondary antibody | 1 vial |
| Mouse uPA standard | 1 vial |
| TMB Substrate Solution | 1 x 10ml |
| uPA Antibody pre-coated 96 well microplate (12 x 8 well strips) | 1 x 96 tests |

関連性

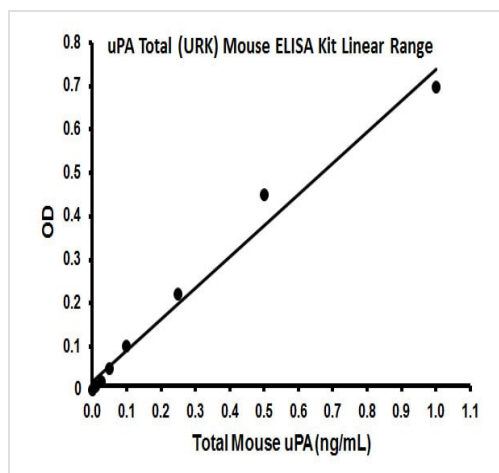
Urokinase type plasminogen activator (uPA) is secreted from cells as a precursor which is activated to the two chain form consisting of an A and B chain. The active high MW form is further processed by removal of an amino terminal fragment to an active low MW form (35 kDa). uPA is a serine protease that activates plasminogen to plasmin. High levels of uPA and plasminogen activator inhibitor type 1 (PAI 1) in breast cancer tissue extracts have been associated with rapid disease progression. The malignant phenotype of prostatic tumor cells correlates with the expression of both uPA and its cell membrane receptor (uPAR).

画像



Full range standard curve using ab198512 uPA Total (URK) Mouse ELISA Kit.

Standard Curve.



Linear range standard curve using ab198512 uPA Total (URK) Mouse 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