

Rat uPA ELISA kit (total uPA antigen) ab272781

画像数 1

医薬用外劇物

製品の概要

製品名	Rat uPA ELISA kit (total uPA antigen)
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Tissue Extracts, Cell culture media, EDTA Plasma, Cit plasma
アッセイタイプ	Quantitative
検出範囲	0.01 ng/ml - 10 ng/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Rat
製品の概要	<p>Rat uPA ELISA kit (total uPA antigen) (ab272781) is intended for the quantitative determination of total urokinase plasminogen activator in rat plasma, serum, urine, cell culture media, or tissue extracts.</p> <p>Rat 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-rat uPA primary antibody binds to the captured enzyme. Excess antibody is washed away and bound polyclonal antibody is then reacted with 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.</p>
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.

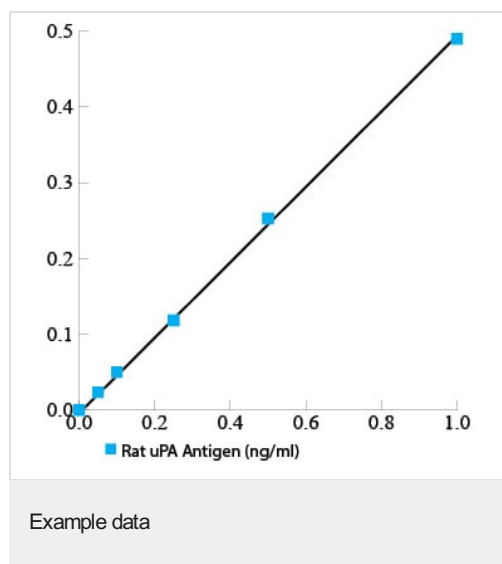
内容	1 x 96 tests
10X Wash Buffer	1 x 50ml
Anti-Rabbit HRP Secondary Reagent	1 vial
Anti-Rat uPA Primary Antibody Lyophilized Vial	1 vial

内容	1 x 96 tests
Rat uPA ELISA Plate	1 unit
Rat uPA Standard Lyophilized Vial	1 vial
TMB Substrate	1 x 10ml

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



A typical standard curve. Example only.

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