

Human Urokinase type Plasminogen Activator ELISA Kit
ab108917

4 References 1 图像

概述

产品名称 人Urokinase type Plasminogen活化剂ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Overall				4.9%

批次间				
样品	n	Mean	SD	CV%
Overall				7.2%

样品类型 Cell culture supernatant, Urine, Serum, Plasma, Tissue

检测类型 Sandwich (quantitative)

灵敏度 = 45 pg/ml

范围 0.125 ng/ml - 4 ng/ml

回收率 98 %

检测时间 4h 00m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Urokinase type Plasminogen Activator Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Urokinase type Plasminogen Activator (uPA) in plasma, serum, urine, milk, cell culture supernatants and tissue extracts.

A Urokinase type Plasminogen Activator specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Urokinase type Plasminogen Activator specific biotinylated detection antibody is added and then followed by washing with wash buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize HRP enzymatic

reaction. TMB is catalyzed by HRP 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 Urokinase type Plasminogen Activator captured in plate.

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

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
70X Biotinylated Human Urokinase type Plasminogen Activator Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml
Urokinase type Plasminogen Activator Microplate (12 x 8 well strips)	1 unit
Urokinase type Plasminogen Activator Standard	1 vial

功能 Specifically cleave the zymogen plasminogen to form the active enzyme plasmin.

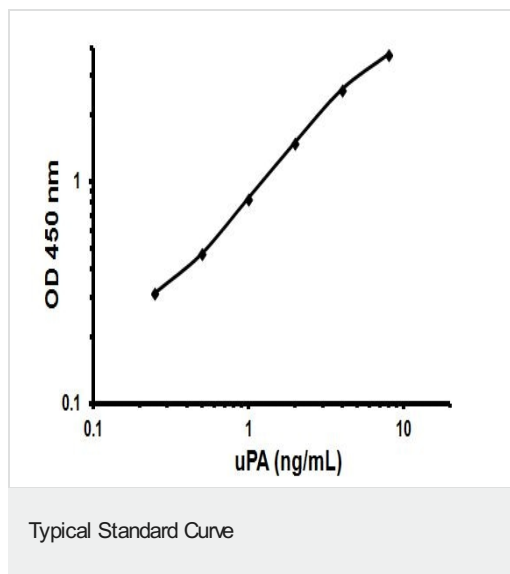
组织特异性 Expressed in the prostate gland and prostate cancers.

序列相似性
Belongs to the peptidase S1 family.
Contains 1 EGF-like domain.
Contains 1 kringle domain.
Contains 1 peptidase S1 domain.

翻译后修饰 Phosphorylation of Ser-158 and Ser-323 abolishes proadhesive ability but does not interfere with receptor binding.

细胞定位 Secreted.

图片



Representative Standard Curve using ab108917.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors