

Human uPA ELISA Kit (URK) ab119611

★★★★★ 1 Abreviews 4 References 1 图像

概述

产品名称	人uPA ELISA试剂盒(URK)				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	1	16	1211pg/ml	8.77	= 4.5%
	2	16	583pg/ml	39.06	= 6.7%
	3	16	1373pg/ml	74.14	= 5.4%
	批次间				
	样品	n	Mean	SD	CV%
	1	24	211pg/ml	12.87	= 6.1%
	2	24	536pg/ml	44.48	= 8.3%
	3	24	1493pg/ml	103.01	= 6.9%
样品类型	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 5 pg/ml				
范围	62.5 pg/ml - 4000 pg/ml				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's Human uPA (URK) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is				

A uPA specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal

antibody from goat specific for uPA is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the 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 Human uPA amount of sample captured in plate.

说明

Plasminogen activator, urokinase (PLAU, uPA) converts plasminogen to plasmin. Plasmin is involved in processing of amyloid precursor protein and degrades secreted and aggregated amyloid-beta, a hallmark of Alzheimer disease (AD). Urokinase has a molecular mass of about 54 kD and is composed of 2 disulfide-linked chains, A and B, of molecular masses 18 kD and 33 kD, respectively. It localized on 10q24. uPA facilitates cell migration by localizing proteolysis on the cell surface and by inducing intracellular signalling pathways. In human vascular smooth muscle cell (VSMC), uPA stimulates migration via the uPA receptor (uPAR) signalling complex containing TYK2 and phosphatidylinositol 3-kinase (PI3-K).

平台

Microplate

性能

存放说明

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

组件	标识符	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human uPA antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human uPA antibody		1 x 100µl
Lyophilized recombinant Human uPA standard		2 x 10ng
Plate Seal		4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

功能

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

组织特异性

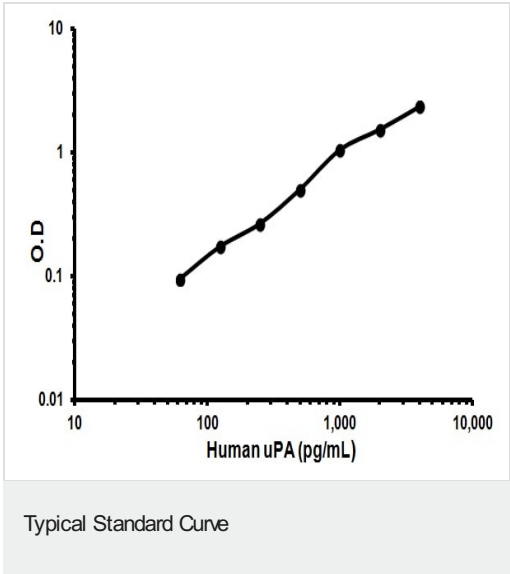
Expressed in the prostate gland and prostate cancers.

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 ab119611

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