

Human PAI1 ELISA Kit ab157528

2 图像

概述

产品名称	人PAI1 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	1	20	2.8U/ml	0.257	9.18%
	2	20	6.82U/ml	0.323	4.74%
	3	20	29.6U/ml	2.57	8.68%
	批次间				
	样品	n	Mean	SD	CV%
	1	10	4.2U/ml	0.399	9.52%
	2	10	8.07U/ml	0.763	9.46%
	3	10	26.8U/ml	2.1	7.85%
样品类型	Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	= 0.11 U/ml				
范围	0.125 U/ml - 100 U/ml				
回收率	96.45 %				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's PAI1 ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for is for the				

PAI1 is involved in the regulation of the blood fibrinolytic system. Increased plasma levels of PAI1 are implicated in the impairment of fibrinolytic function and may be associated with thrombotic

diseases. Levels of PAI1 increase with age and are elevated in conditions such as normal pregnancy and sepsis.

Functionally active PAI1 present in plasma reacts with urokinase coated and dried on a microtiter plate. Latent or complexed PAI1 will not bind to the plate or be detected. Unbound PAI1 samples are aspirated and an anti-PAI1 primary antibody is added. Excess primary antibody is then aspirated. The bound antibody, which is proportional to the original active PAI1 present in the samples, is then reacted with the horseradish peroxidase conjugated secondary antibody. Following an additional washing step, TMB substrate solution is then used for color development at 450nm. The amount of color development is directly proportional to the concentration of active PAI1 in the sample.

Get results in 90 minutes with Human PAI1 ELISA Kit (SERPINE1) ([ab184863](#)) from our SimpleStep ELISA® range.

平台 Microplate

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Wash Buffer	1 x 50ml
Anti-Human PAI1 Primary Antibody	1 vial
Assay Diluent	1 x 10ml
HRP secondary antibody	1 vial
Human PAI1 high activity standard (260U)	1 vial
Human PAI1 zero unit activity standard (0 U)	2 vials
TMB Substrate Solution	1 x 10ml
uPA Antibody pre-coated 96 well microplate (12 x 8 well strips)	1 unit

功能 This inhibitor acts as 'bait' for tissue plasminogen activator, urokinase, and protein C. Its rapid interaction with TPA may function as a major control point in the regulation of fibrinolysis.

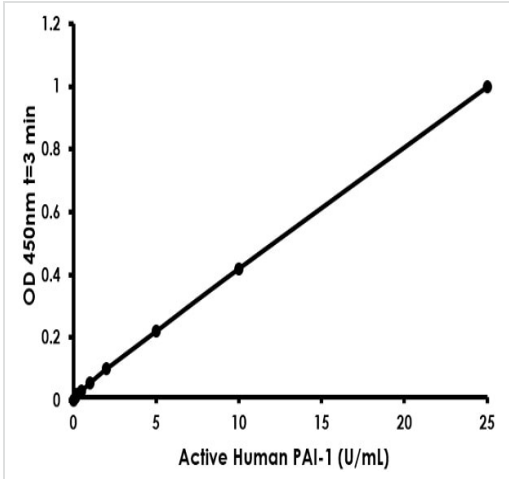
组织特异性 Found in plasma and platelets and in endothelial, hepatoma and fibrosarcoma cells.

疾病相关 Defects in SERPINE1 are the cause of plasminogen activator inhibitor-1 deficiency (PAI-1D) [MIM:613329]. It is a hematologic disorder characterized by increased bleeding after trauma, injury, or surgery. Affected females have menorrhagia. The bleeding defect is due to increased fibrinolysis of fibrin blood clots due to deficiency of plasminogen activator inhibitor-1, which inhibits tissue and urinary activators of plasminogen.
Note=High concentrations of SERPINE1 seem to contribute to the development of venous but not arterial occlusions.

序列相似性 Belongs to the serpin family.

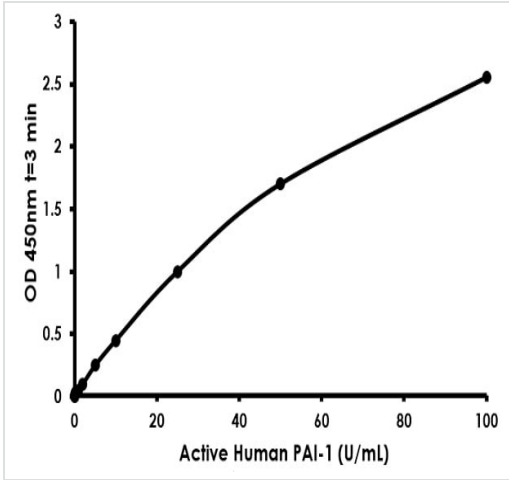
翻译后修饰	Inactivated by proteolytic attack of the urokinase-type (u-PA) and the tissue-type (TPA), cleaving the 369-Arg-Met-370 bond.
细胞定位	Secreted.

图片



Typical Standard Curve for ab157528 PAI-1 Human ELISA Linear Range

Typical Standard Curve for ab157528 PAI-1 Human ELISA Linear Range



Typical Standard Curve for ab157528 PAI-1 Human ELISA Full range

Typical Standard Curve for ab157528 PAI-1 Human ELISA Full range

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