

3 References 画像数 2

医薬用外劇物

製品の概要

製品名	Rat PAI1 ELISA Kit (SERPINE1)
-----	-------------------------------

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	0.957ng/ml	0.088	9.24%
Known Conc	20	2.25ng/ml	0.134	5.97%
Known Conc	20	15.6ng/ml	1.01	6.48%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Known Conc	10	0.926ng/ml	0.113	12.2%
Known Conc	10	1.4ng/ml	0.096	6.86%
Known Conc	10	9.17ng/ml	0.858	9.36%

サンプルの種類 Serum, Plasma

アッセイタイプ Quantitative

検出感度 0.079 ng/ml

検出範囲 0.1 ng/ml - 50 ng/ml

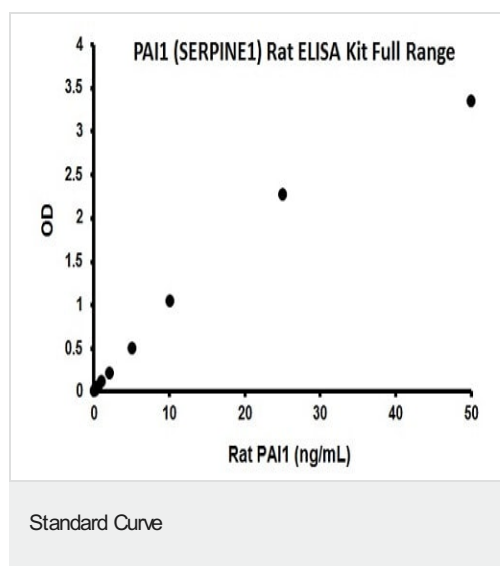
添加回收試驗	104 %
--------	-------

特定サンプルでの回収試験

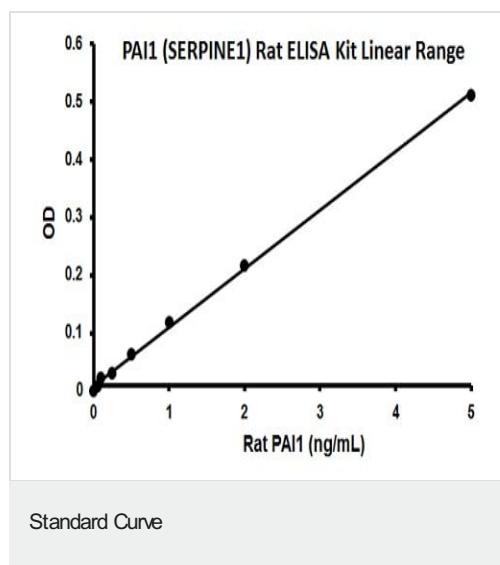
サンプルの種類	平均 %	測定範囲
Spike	104	97% - 108%

ステップ Multiple steps standard assay

種交差性	交差種: Rat 非交差種: Sheep
製品の概要	<p>Abcam's PAI1 (SERPINE1) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of active PAI1 (SERPINE1) in rat plasma and serum.</p> <p>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 and will not be detected. Unbound PAI1 samples are washed away and an anti-PAI1 primary antibody is added. Excess primary antibody is washed away and bound antibody, which is proportional to the original active PAI1 present in the samples, is then bound by the horseradish peroxidase secondary antibody. Following an additional washing step, TMB 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.</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 horseradish peroxidase secondary antibody	1 vial
Anti-rat PAI-1 primary antibody	1 vial
Rat PAI-1 activity 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
機能	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.
組織特異性	Found in plasma and platelets and in endothelial, hepatoma and fibrosarcoma cells.
関連疾患	<p>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.</p> <p>Note=High concentrations of SERPINE1 seem to contribute to the development of venous but not arterial occlusions.</p>
配列類似性	Belongs to the serpin family.
翻訳後修飾	Inactivated by proteolytic attack of the urokinase-type (u-PA) and the tissue-type (TPA), cleaving the 369-Arg-Met-370 bond.
細胞内局在	Secreted.



Full range standard curve using ab198509 PAI1 (SERPINE1) Rat ELISA Kit.



Linear range standard curve using ab198509 PAI1 (SERPINE1) Rat ELISA Kit.

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