

Human Antithrombin III ELISA Kit ab108801

★★★★★ 1 Abreviews 2 References 画像数 7

製品の概要

製品名	Human Antithrombin III ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				2.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.4%
サンプルの種類	Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 0.71 ng/ml			
検出範囲	6.3 ng/ml - 82.5 ng/ml			
添加回収試験	94 %			
全工程の試験時間	4h 0m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's Antithrombin III Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Antithrombin III in urine, milk, saliva, cerebrospinal fluid, plasma, and serum.			

An antithrombin III specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an antithrombin III specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase 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 antithrombin III captured in plate.

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

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	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
40X Biotinylated Human Antithrombin III Antibody	1 x 150µl
Antithrombin III Microplate (12 x 8 well strips)	1 unit
Antithrombin III Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Most important serine protease inhibitor in plasma that regulates the blood coagulation cascade. AT-III inhibits thrombin as well as factors IXa, Xa and XIa. Its inhibitory activity is greatly enhanced in the presence of heparin.

組織特異性

Found in plasma.

関連疾患

Defects in SERPINC1 are the cause of antithrombin III deficiency (AT3D) [MIM:613118]. AT3D is an important risk factor for hereditary thrombophilia, a hemostatic disorder characterized by a tendency to recurrent thrombosis. AT3D is classified into 4 types. Type I: characterized by a 50% decrease in antigenic and functional levels. Type II: has defects affecting the thrombin-binding domain. Type III: alteration of the heparin-binding domain. Plasma AT-III antigen levels are normal in type II and III. Type IV: consists of miscellaneous group of unclassifiable mutations.

配列類似性

Belongs to the serpin family.

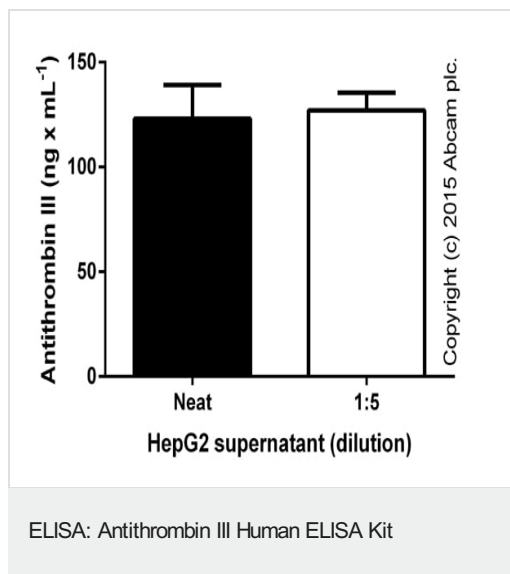
翻訳後修飾

Phosphorylation sites are present in the extracellular medium.

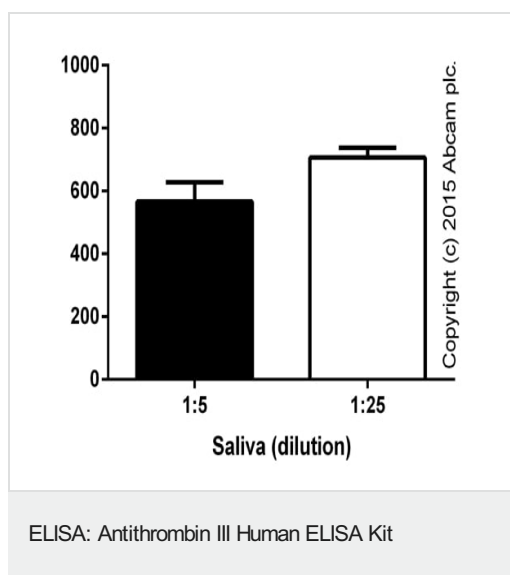
細胞内局在

Secreted > extracellular space.

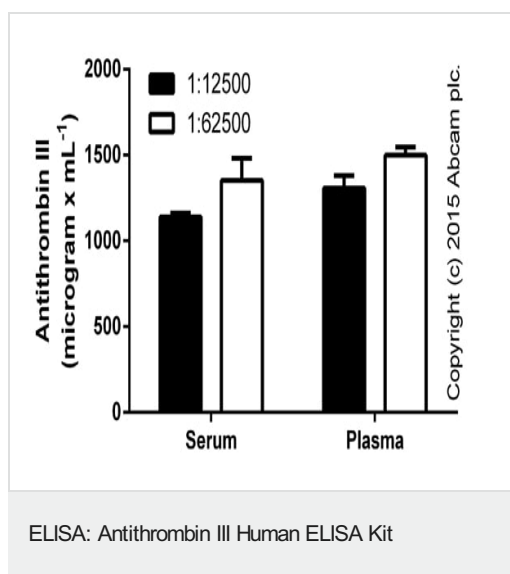
画像



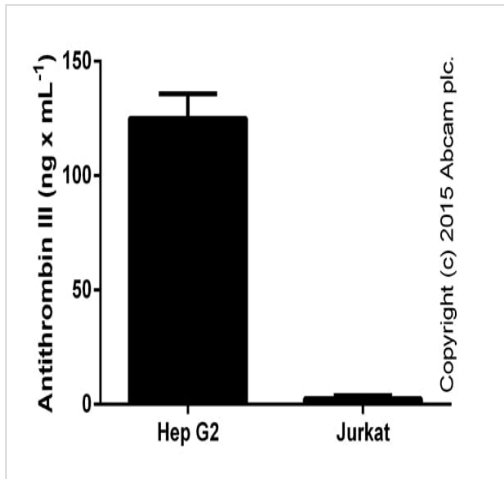
Linearity: Antithrombin III in samples at different dilutions, corrected for the factor of dilution (duplicates +/- SD).



Linearity: Antithrombin III in samples at different dilutions, corrected for the factor of dilution (duplicates +/- SD), showing ng x mL⁻¹ on y-axis.

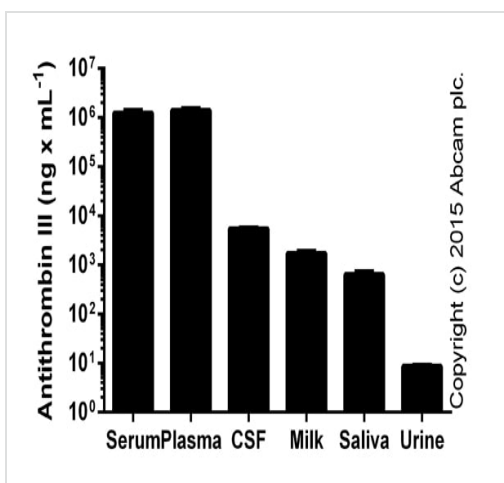


Linearity: Antithrombin III in samples at different dilutions, corrected for the factor of dilution (duplicates +/- SD).



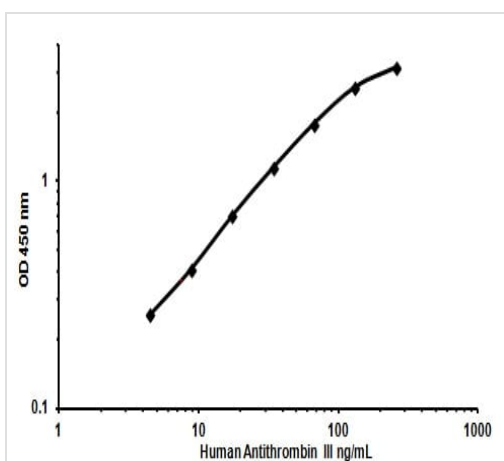
ELISA: Antithrombin III Human ELISA Kit

Antithrombin III in cell culture supernatants (duplicates +/- SD), tested in the dilution range of 1:1-1:100.



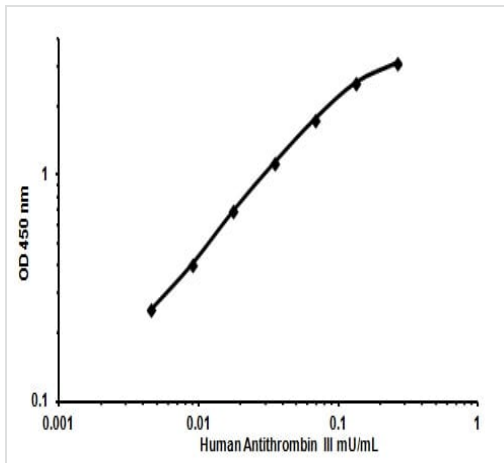
ELISA: Antithrombin III Human ELISA Kit

Antithrombin III in human biological fluids (duplicates +/- SD). The samples were differently diluted (serum and Li-heparine plasma: 1e4-1e5; CSF: 1e2-1e3; milk: 10-100; saliva: 5-100; urine: neat).



Typical Standard Curve

Representative standard curve using ab108801.



Representative standard curve using ab108801.

Typical Standard Curve

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