

Human Thrombin-Antithrombin Complex ELISA Kit (TAT)
ab108907

★★★★☆ 3 Abreviews 18 References 画像数 2

製品の概要

製品名	Human Thrombin-Antithrombin Complex ELISA Kit (TAT)			
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%
				5.4%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				10.4%
サンプルの種類	Cell culture supernatant, Milk, Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 0.78 ng/ml			
検出範囲	1.48 ng/ml - 120 ng/ml			
添加回収試験	96 %			
全工程の試験時間	4h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Thrombin-Antithrombin Complex (TAT) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Thrombin-Antithrombin Complex concentrations in plasma, milk and cell culture supernatants.			

A Thrombin-Antithrombin Complex specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Thrombin-Antithrombin Complex specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate 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 Thrombin-Antithrombin Complex captured in plate.

Note: It is not recommended to perform spike recovery **in plasma samples** due to large amount of endogenous Anti-thrombin III (mg/ml range) which will saturate the capture antibody and interfere with the binding of the TAT standard

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
50X Biotinylated Human Thrombin-Antithrombin Complex Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml
Thrombin-Antithrombin Complex Microplate (12 x 8 well strips)	1 unit
Thrombin-Antithrombin Complex Standard	1 vial

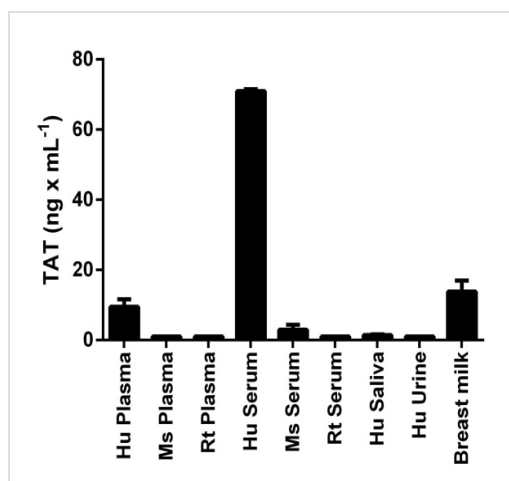
関連性

Thrombin, which cleaves bonds after Arg and Lys, converts fibrinogen to fibrin and activates factors V, VII, VIII, XIII, and, in complex with thrombomodulin, protein C. Functions in blood homeostasis, inflammation and wound healing. Antithrombin: Most important serine protease inhibitor in plasma that regulates the blood coagulation cascade. Antithrombin inhibits thrombin, matriptase-3/TMPRSS7, as well as factors IXa, Xa and XIa. Its inhibitory activity is greatly enhanced in the presence of heparin.

細胞内局在

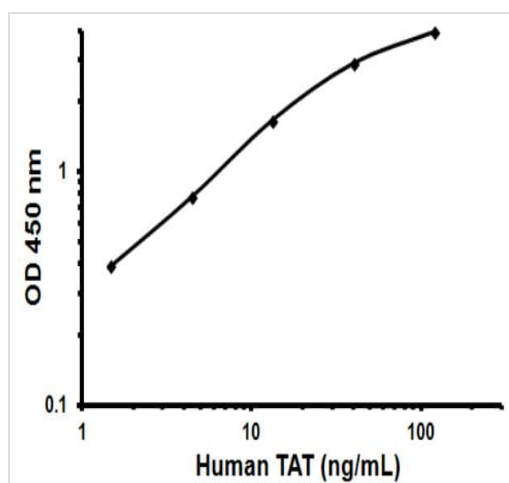
Secreted

画像



TAT measured in biological fluids, with background signal subtracted (duplicates +/- SD).

Sandwich ELISA - Thrombin-Antithrombin Complex
(TAT) Human ELISA Kit



Representative Standard Curve Using ab108907.

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