

Mouse Antithrombin III ELISA Kit (SERPINC1) ab108800

1 References 2 图像

概述

产品名称	小鼠Antithrombin III ELISA试剂盒(SERPINC1)			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Overall			6.2%
	批次间			
	样品	n	Mean	SD
	Overall			10.7%
样品类型	Cell culture supernatant, Urine, Serum, Plasma			
检测类型	Sandwich (quantitative)			
灵敏度	= 1.7 ng/ml			
范围	6.25 ng/ml - 50 ng/ml			
回收率	96 %			
检测时间	4h 00m			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Mouse			
产品概述	Abcam's Antithrombin III (SERPINC1) Mouse <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Antithrombin III concentrations in mouse plasma, serum, urine and cell culture supernatants.			

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.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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 Mouse Antithrombin III Antibody	1 x 120µl
Antithrombin III Standard	1 vial
Antithrombin III Microplate (12 x 8 well strips)	1 unit
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

功能 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.

组织特异性 Found in plasma.

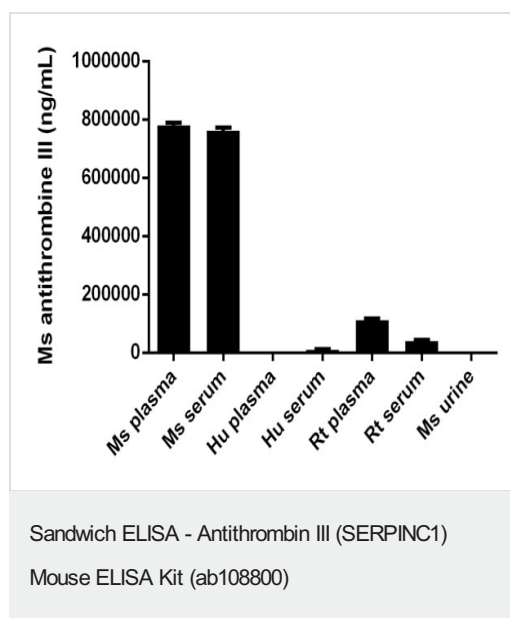
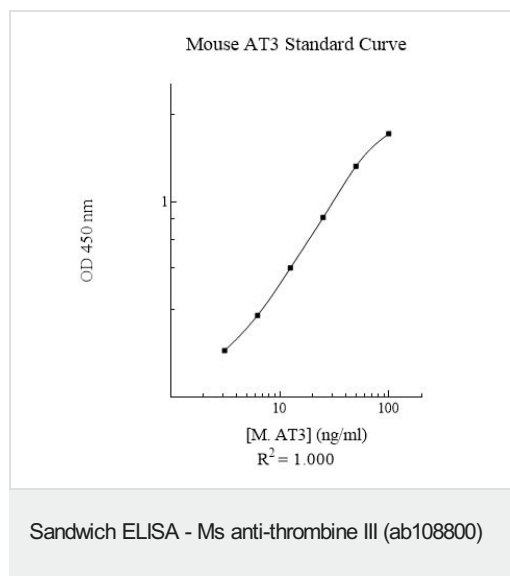
疾病相关 Defects in SERPINC1 are the cause of antithrombin III deficiency (AT3D) [MIM:613118]. AT3D is an important risk factor for hereditary thrombophilia, a hemostatic disorder characterized by a tendency to recurrent thrombosis. AT3D is classified into 4 types. Type I: characterized by a 50% decrease 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.

序列相似性 Belongs to the serpin family.

翻译后修饰 Phosphorylation sites are present in the extracellular medium.

细胞定位 Secreted > extracellular space.

图片



Ms antithrombine III measured in biological fluids showing quantity (ng) per mL of sample tested

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