

Mouse Antithrombin III ELISA Kit (SERPINC1) ab108800

1 References 2 Images

Overview

Product name	Mouse Antithrombin III ELISA Kit (SERPINC1)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				6.2%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				10.7%
Sample type	Cell culture supernatant, Urine, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	= 1.7 ng/ml				
Range	6.25 ng/ml - 50 ng/ml				
Recovery	96 %				
Assay time	4h 00m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	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.

Platform Microplate

Properties

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

Components	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

Function 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.

Tissue specificity Found in plasma.

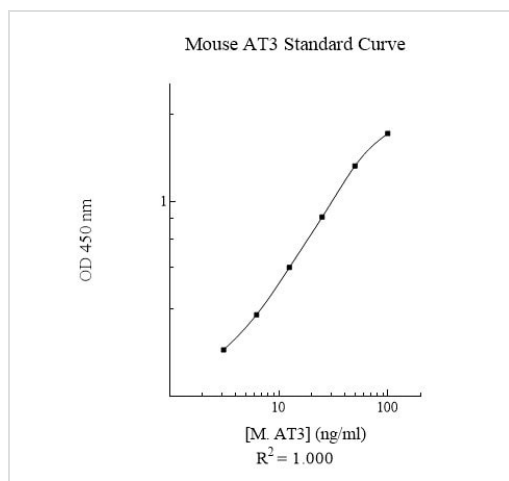
Involvement in disease Defects in SERPINC1 are the cause of antithrombin III deficiency (AT3D) [MIM:613118]. AT3D is an important risk factor for hereditary thrombophilia, a hemostatic disorder characterized by a tendency to recurrent thrombosis. AT3D is classified into 4 types. Type I: characterized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

Sequence similarities Belongs to the serpin family.

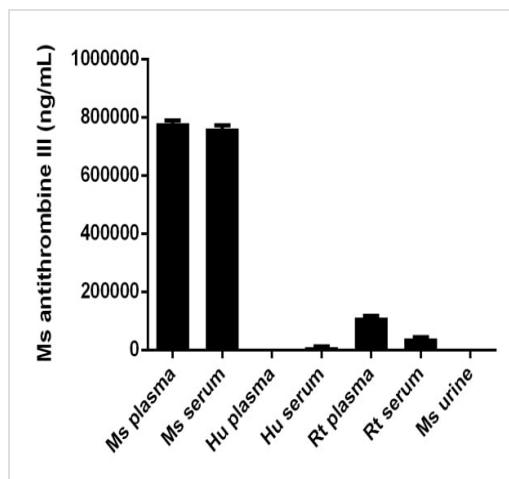
Post-translational modifications Phosphorylation sites are present in the extracellular medium.

Cellular localization Secreted > extracellular space.

Images



Sandwich ELISA - Ms anti-thrombine III (ab108800)



Sandwich ELISA - Antithrombin III (SERPINC1)
Mouse ELISA Kit (ab108800)

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors