

Human Protein S ELISA Kit ab125969

1 References 1 图像

概述

产品名称	人蛋白S ELISA试剂盒			
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			9%
样品类型	Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate			
检测类型	Competitive			
灵敏度	0.22 µg/ml			
范围	1 µg/ml - 4 µg/ml			
回收率	97 %			
检测时间	4h 00m			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	Abcam's Protein S Human <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Human protein S in plasma, serum, cell culture supernatants cell lysate and tissue samples.			

A Protein S specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Protein S 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 inversely proportional to the amount of Protein S 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
1X Biotinylated Human Protein S (Lyophilized)	1 vial
20X Wash Buffer Concentrate	1 x 30ml
Chromogen Substrate	1 x 7ml
Protein S Microplate (12 x 8 well strips)	1 unit
Protein S Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

功能 Anticoagulant plasma protein; it is a cofactor to activated protein C in the degradation of coagulation factors Va and VIIIa. It helps to prevent coagulation and stimulating fibrinolysis.

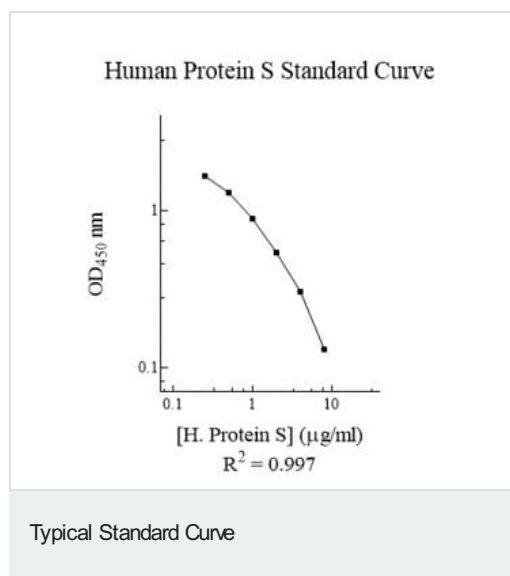
组织特异性 Plasma.

疾病相关 Defects in PROS1 are the cause of protein S deficiency (PROS1D) [MIM:612336]; also known as thrombophilia due to protein S deficiency. PROS1D is a cause of hereditary thrombophilia, a hemostatic disorder characterized by impaired regulation of blood coagulation and a tendency to recurrent venous thrombosis. However, many adults with heterozygous disease may be asymptomatic. Based on the plasma levels of total and free PROS1 antigen as well as the serine protease-activated protein C cofactor activity, three types of PROS1D have been described: type I, characterized by reduced total and free PROS1 antigen levels together with reduced anticoagulant activity; type III, in which only free PROS1 antigen and PROS1 activity levels are reduced; and the rare type II which is characterized by normal concentrations of both total and free PROS1 antigen, but low cofactor activity.

序列相似性 Contains 4 EGF-like domains.
Contains 1 Gla (gamma-carboxy-glutamate) domain.
Contains 2 laminin G-like domains.

翻译后修饰 The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.

细胞定位 Secreted.



Representative Standard Curve using ab125969

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