

Human Thrombomodulin ELISA Kit (CD141) ab46508

★★★★★ [1 Abreviews](#) [18 References](#) [画像数 2](#)

製品の概要

製品名	Human Thrombomodulin ELISA Kit (CD141)			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Serum	6		CV%
				3.9%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Serum	6		CV%
				9.8%
サンプルの種類	Cell culture supernatant, Serum, Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 0.31 ng/ml			
検出範囲	0.625 ng/ml - 20 ng/ml			
添加回収試験	109 %			
全工程の試験時間	1h 45m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's Thrombomodulin (CD141) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Thrombomodulin (CD141) in serum, plasma, buffered solutions and supernatants.			

A monoclonal antibody specific for Thrombomodulin has been coated onto the wells of the microtiter strips provided. Samples, including standards of known Thrombomodulin concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for Thrombomodulin are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of Thrombomodulin present in the samples.

This kit will recognize both endogenous and recombinant Human Thrombomodulin.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	1 x 96 tests	2 x 96 tests	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml	1 x 7ml	1 x 13ml
Biotinylated anti-Thrombomodulin	Red	1 x 400µl	2 x 400µl	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml	1 x 11ml	1 x 24ml
Control	Silver	2 vials	4 vials	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml	1 x 12ml	1 x 23ml
Stop Reagent	Black	1 x 11ml	2 x 11ml	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl	2 x 5µl	4 x 5µl
Thrombomodulin Microplate (12 x 8 well strips)		1 unit	2 units	1 unit	2 units
Thrombomodulin Standard (lyophilized)	Yellow	2 vials	4 vials	2 vials	4 vials

機能

Thrombomodulin is a specific endothelial cell receptor that forms a 1:1 stoichiometric complex with thrombin. This complex is responsible for the conversion of protein C to the activated protein C (protein Ca). Once evolved, protein Ca scissions the activated cofactors of the coagulation mechanism, factor Va and factor VIIIa, and thereby reduces the amount of thrombin generated.

組織特異性

Endothelial cells are unique in synthesizing thrombomodulin.

関連疾患

Defects in THBD are the cause of thrombophilia due to thrombomodulin defect (THR-THBD) [MIM:188040]. A hemostatic disorder characterized by a tendency to thrombosis.

Defects in THBD are a cause of susceptibility to hemolytic uremic syndrome atypical type 6 (AHUS6) [MIM:612926]. An atypical form of hemolytic uremic syndrome. It is a complex genetic disease characterized by microangiopathic hemolytic anemia, thrombocytopenia, renal failure and absence of episodes of enterocolitis and diarrhea. In contrast to typical hemolytic uremic syndrome, atypical forms have a poorer prognosis, with higher death rates and frequent progression to end-stage renal disease. Note=Susceptibility to the development of atypical hemolytic uremic syndrome can be conferred by mutations in various components of or regulatory factors in the complement cascade system. Other genes may play a role in modifying the phenotype.

配列類似性

Contains 1 C-type lectin domain.

Contains 6 EGF-like domains.

翻訳後修飾

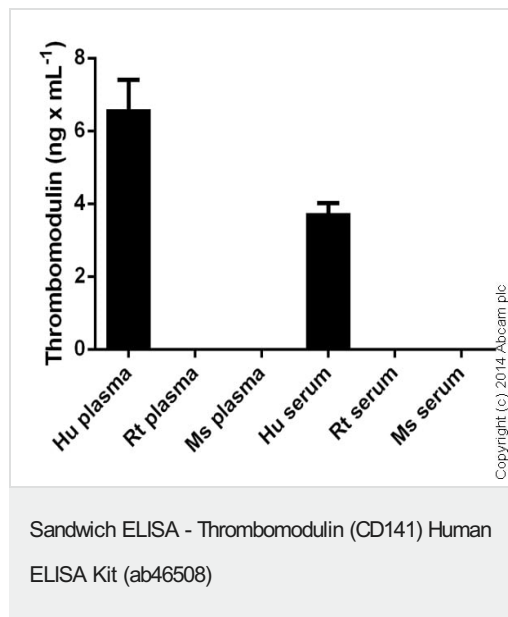
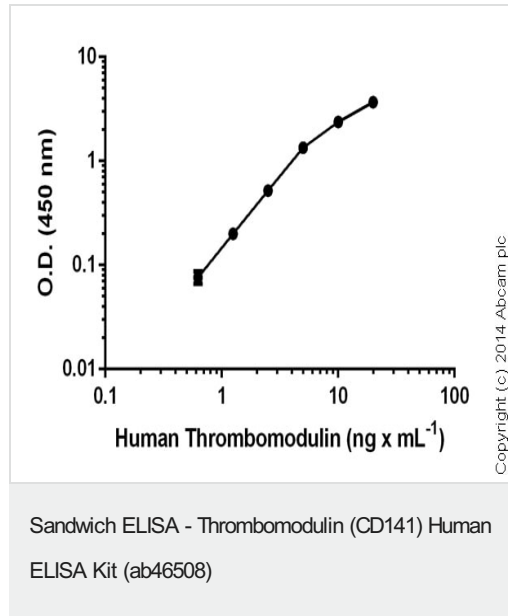
N-glycosylated.

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

細胞内局在

Membrane.

画像



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