

Human ADAMTS1 ELISA Kit ab213751

1 图像

概述

产品名称

人ADAMTS1 ELISA试剂盒

检测方法

Colorimetric

精确度

批次内

样品	n	Mean	SD	CV%
1	16	145pg/ml	8.84	= 6.1%
2	16	681pg/ml	37.45	= 5.5%
3	16	2886pg/ml	207.79	= 7.2%

批次间

样品	n	Mean	SD	CV%
1	24	132pg/ml	10.56	= 8%
2	24	732pg/ml	51.97	= 7.1%
3	24	2843pg/ml	213.22	= 7.5%

样品类型

Cell culture supernatant, Serum, Hep Plasma

检测类型

Sandwich (quantitative)

灵敏度

< 10 pg/ml

范围

93.8 pg/ml - 6000 pg/ml

检测时间

3h 30m

实验步骤

Multiple steps standard assay

种属反应性

与反应: Human

产品概述

The Human ADAMTS1 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213751) is designed for the quantitative measurement of Human ADAMTS1 in cell culture supernatants, serum and plasma (heparin).

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from mouse specific for ADAMTS1 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: F253-A734) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for ADAMTS1 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human ADAMTS1 amount of sample captured in plate.

说明 ADAMTS1 (A Disintegrin-Like and Metalloproteinase with Thrombospondin Type 1 Motif, 1), also known as METH1, is an enzyme that in humans is encoded by the ADAMTS1 gene. ADAMTS is a family of proteins believed to be anchored to the extracellular matrix (ECM) through interactions with aggrecan or other matrix components by one or more thrombospondin type 1 motifs. The ADAMTS1 gene is mapped to 21q21.2 based on sequence similarity between the ADAMTS1 sequence and a chromosome 21q21.2 clone. It has been found that ADAMTS1 can disrupt angiogenesis *in vivo* and *in vitro* more efficiently than ADAMTS8, THBS1, or endostatin.

平台 Pre-coated microplate (12 x 8 well strips)

性能

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

组件	标识符	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human ADAMTS1 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Human ADAMTS1 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human ADAMTS1 standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

功能 Cleaves aggrecan, a cartilage proteoglycan, and may be involved in its turnover (By similarity). Has angiogenic inhibitor activity. Active metalloprotease, which may be associated with various inflammatory processes as well as development of cancer cachexia. May play a critical role in follicular rupture.

序列相似性 Contains 1 disintegrin domain.

Contains 1 peptidase M12B domain.

Contains 3 TSP type-1 domains.

结构域

The spacer domain and the TSP type-1 domains are important for a tight interaction with the extracellular matrix.

The conserved cysteine present in the cysteine-switch motif binds the catalytic zinc ion, thus inhibiting the enzyme. The dissociation of the cysteine from the zinc ion upon the activation-peptide release activates the enzyme.

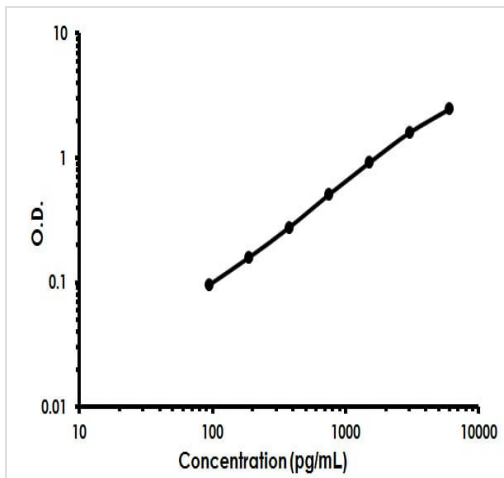
翻译后修饰

The precursor is cleaved by a furin endopeptidase.

细胞定位

Secreted > extracellular space > extracellular matrix.

图片



Human ADAMTS1 ELISA Kit (ab213751) Standard Curve.

Human ADAMTS1 ELISA Kit (ab213751) 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.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