

Rat TIMP-2 ELISA Kit ab213923

★★★★★ 1 Abreviews 1 图像

概述

产品名称	大鼠TIMP-2 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	1	16	240pg/ml	10.56	4.4%
	2	16	1342pg/ml	68.44	9.3%
	3	16	4533pg/ml	339.97	7.5%
	批次间				
	样品	n	Mean	SD	CV%
	1	24	227pg/ml	10.21	4.5%
	2	24	1463pg/ml	95.09	6.5%
	3	24	4769pg/ml	443.51	9.3%
样品类型	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate				
检测类型	Sandwich (quantitative)				
灵敏度	< 10 pg/ml				
范围	156 pg/ml - 10000 pg/ml				
检测时间	3h 30m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Rat				
产品概述	The Rat TIMP-2 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213923) is designed for				

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A polyclonal antibody from goat specific for TIMP-2 has been precoated onto 96-well plates. Standards and test samples are added to the wells; a biotinylated detection polyclonal antibody from goat specific for TIMP-2 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was 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 Rat TIMP-2 amount of sample captured in plate.

**说明** TIMP-2 gene is encoded by 5 exons spanning 83 kb of genomic DNA. TIMP-2 is 83 kilobase pairs (kb) long with exon-intron splicing sites located in preserved positions among the three members of the TIMP family. The gene for tissue inhibitor of metalloproteinases-2 is localized on human chromosome arm 17q25. TIMP-2 abrogates angiogenic factor-induced endothelial cell proliferation *in vitro* and angiogenesis *in vivo* independent of MMP inhibition.

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

**功能** Complexes with metalloproteinases (such as collagenases) and irreversibly inactivates them by binding to their catalytic zinc cofactor. Known to act on MMP-1, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, MMP-10, MMP-13, MMP-14, MMP-15, MMP-16 and MMP-19.

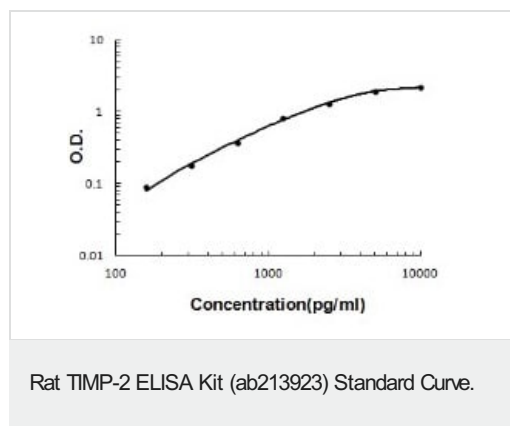
**序列相似性** Belongs to the protease inhibitor I35 (TIMP) family.  
Contains 1 NTR domain.

**翻译后修饰** The activity of TIMP2 is dependent on the presence of disulfide bonds.

**细胞定位** Secreted.

## 图片

---



Rat TIMP-2 ELISA Kit (ab213923) 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