

Mouse TIMP1 ELISA Kit ab100745

2 References 画像数 1

製品の概要

製品名	Mouse TIMP1 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture extracts, Tissue Extracts
アッセイタイプ	Sandwich (quantitative)
検出感度	< 30 pg/ml
検出範囲	24.69 pg/ml - 18000 pg/ml
添加回収試験	100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture extracts	114	105% - 123%
Tissue Extracts	106.5	94% - 116%

ステップ	Multiple steps standard assay
種交差性	交差種: Mouse

**製品の概要**

Abcam's TIMP1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme linked immunosorbent assay for the quantitative measurement of Mouse TIMP1 in cell lysates and tissue lysates.

This assay employs an antibody specific for Mouse TIMP1 coated on a 96-well plate. Standards and samples are pipetted into the wells and TIMP1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse TIMP1 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of TIMP1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**特記事項**

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム	Microplate
------------	------------

## 製品の特性

### 保存方法

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

内容	1 x 96 tests
120X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
2X Cell Lysis Buffer	1 x 5ml
5X Assay Diluent	1 x 15ml
5X Sample Diluent Buffer	1 x 10ml
Biotinylated anti-Mouse TIMP1	2 vials
Recombinant Mouse TIMP1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TIMP1 Microplate (12 x 8 wells)	1 unit
TMB One-Step Substrate Reagent	1 x 12ml

### 機能

Complexes with metalloproteinases (such as collagenases) and irreversibly inactivates them by binding to their catalytic zinc cofactor. Also mediates erythropoiesis in vitro; but, unlike IL-3, it is species-specific, stimulating the growth and differentiation of only human and murine erythroid progenitors. Known to act on MMP-1, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, MMP-10, MMP-11, MMP-12, MMP-13 and MMP-16. Does not act on MMP-14.

### 配列類似性

Belongs to the protease inhibitor I35 (TIMP) family.  
Contains 1 NTR domain.

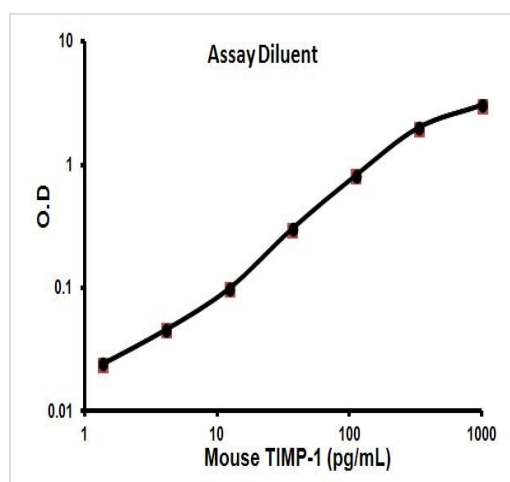
### 翻訳後修飾

The activity of TIMP1 is dependent on the presence of disulfide bonds.

### 細胞内局在

Secreted.

## 画像



Representative Standard Curve using ab100745

Representative 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.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