

### Mouse MMP3 ELISA Kit ab100731

2 References [画像数 1](#)

#### 製品の概要

製品名	Mouse MMP3 ELISA Kit
検出方法	Colorimetric
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	< 10 pg/ml
検出範囲	13.72 pg/ml - 10000 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 supernatant	117.4	106% - 130%
Serum	120.6	101% - 140%
Plasma	111.6	101% - 137%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Abcam's MMP3 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse MMP3 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse MMP3 coated on a 96-well plate. Standards and samples are pipetted into the wells and MMP3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse MMP3 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 MMP3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated anti-Mouse MMP-3	2 vials
MMP-3 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse MMP-3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**機能**

Can degrade fibronectin, laminin, gelatins of type I, III, IV, and V; collagens III, IV, X, and IX, and cartilage proteoglycans. Activates procollagenase.

**配列類似性**

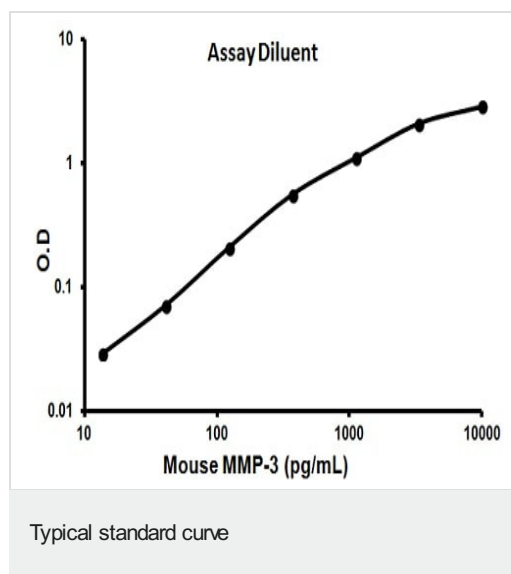
Belongs to the peptidase M10A family.  
Contains 4 hemopexin-like domains.

**ドメイン**

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.

**細胞内局在**

Secreted > extracellular space > extracellular matrix.

**画像**

Representative standard curve using ab100731

**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