

Human MMP10 ELISA Kit ab100602

1 图像

概述

产品名称	人MMP10 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 3 pg/ml
范围	2.06 pg/ml - 1500 pg/ml
回收率	96 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	98.31	87% - 109%
Serum	96.54	85% - 108%
Plasma	94.57	84% - 105%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's MMP10 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human MMP10 pro and active forms in serum, plasma (Collect plasma using heparin as an anticoagulant. EDTA and Citrate are not recommended), and cell culture supernates.

This assay employs an antibody specific for Human MMP10 coated on a 96- well plate. Standards and samples are pipetted into the wells and MMP10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human MMP10 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 MMP10 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明 Optimisation may be required with urine samples.

平台 Microplate

性能

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200μl
Biotinylated anti-Human MMP10	2 vials
MMP10 Microplate (12 x 8 wells)	1 unit
Recombinant Human MMP10 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

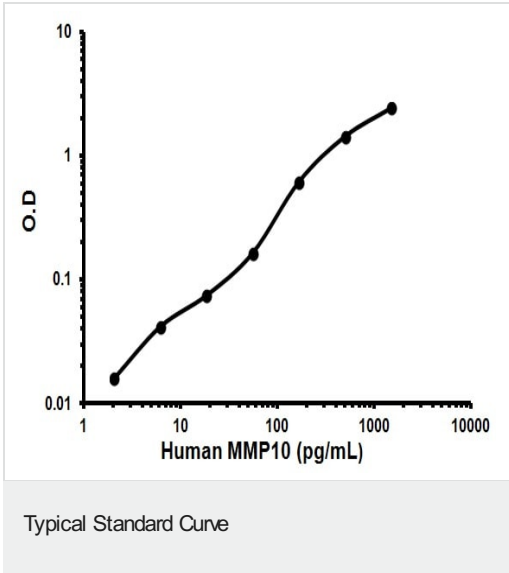
功能Can degrade fibronectin, gelatins of type I, III, IV, and V; weakly collagens III, IV, and V. Activates procollagenase.

序列相似性Belongs to the peptidase M10A family.  
Contains 4 hemopexin-like domains.

结构域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.

细胞定位Secreted > extracellular space > extracellular matrix.

图片



Representative Standard Curve using ab100602.

**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