

Human MMP10 ELISA Kit ab100602

画像数 1

製品の概要

製品名	Human MMP10 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	< 3 pg/ml
検出範囲	2.06 pg/ml - 1500 pg/ml
添加回収試験	96 %

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	98.31	87% - 109%
Serum	96.54	85% - 108%
Plasma	94.57	84% - 105%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

特記事項 Optimisation may be required with urine samples.

試験プラットフォーム Microplate

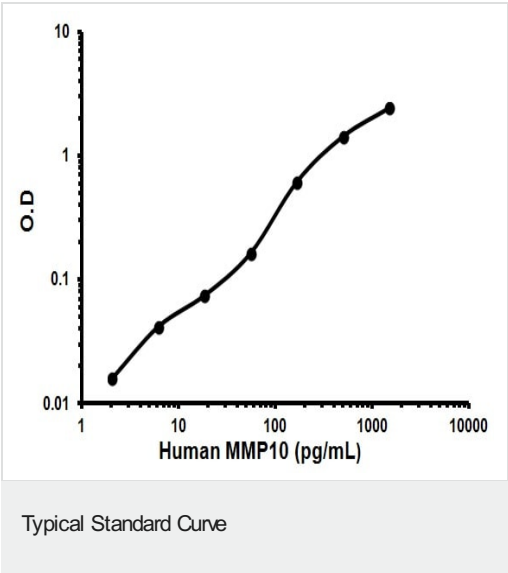
製品の特性

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
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

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