

Human MMP10 ELISA Kit ab100602

1 Image

Overview

|                  |   |
|------------------|---|
| Product name     | Human MMP10 ELISA Kit                   |
| Detection method | Colorimetric                            |
| Sample type      | Cell culture supernatant, Serum, Plasma |
| Assay type       | Sandwich (quantitative)                 |
| Sensitivity      | < 3 pg/ml                               |
| Range            | 2.06 pg/ml - 1500 pg/ml                 |
| Recovery         | 96 %                                    |

Sample specific recovery

| Sample type              | Average % | Range      |
|--------------------------|-----------|------------|
| Cell culture supernatant | 98.31     | 87% - 109% |
| Serum                    | 96.54     | 85% - 108% |
| Plasma                   | 94.57     | 84% - 105% |

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview 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.

Notes Optimisation may be required with urine samples.

Platform Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

| Components                                     | 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     |

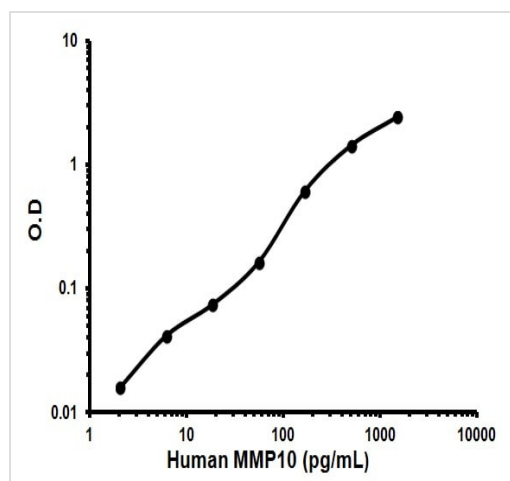
**Function** Can degrade fibronectin, gelatins of type I, III, IV, and V; weakly collagens III, IV, and V. Activates procollagenase.

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

**Domain** 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.

**Cellular localization** Secreted > extracellular space > extracellular matrix.

## Images



Representative Standard Curve using ab100602.

Typical 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.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors