

Human MMP3 ELISA Kit ab100607

★★★★★ 1 Abreviews 4 References 1 Image

Overview

Product name	Human MMP3 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.3 ng/ml
Range	0.41 ng/ml - 100 ng/ml
Recovery	95 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	94.72	84% - 103%
Serum	95.39	83% - 103%
Plasma	96.25	85% - 104%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Abcam’s MMP3 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human MMP3 pro and active forms in serum, plasma, and cell culture supernatants. (Collect plasma using heparin as an anticoagulant. EDTA and Citrate are not recommended).

This assay employs an antibody specific for Human 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-Human 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.

Get results in 90 minutes with Human MMP3 ELISA Kit (ab189572) from our SimpleStep ELISA®

range.

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
1400X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human MMP3	2 vials
MMP3 Microplate (12 strips x 8 wells)	1 unit
MMP3 Standard Human (recombinant) (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

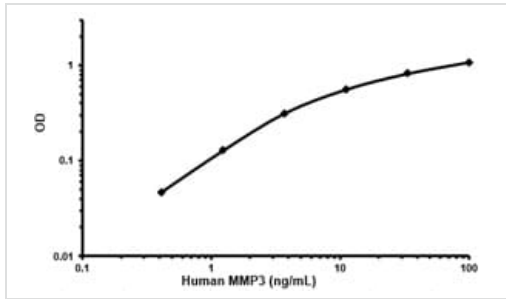
Function Can degrade fibronectin, laminin, gelatins of type I, III, IV, and V; collagens III, IV, X, and IX, and cartilage proteoglycans. Activates procollagenase.

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 ab100607.

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