

Product datasheet

Human MMP1 Matched Antibody Pair Kit ab218184

1 Image

Overview

<b>Product name</b>	Human MMP1 Matched Antibody Pair Kit
<b>Detection method</b>	Colorimetric
<b>Assay type</b>	ELISA set
<b>Range</b>	125 pg/ml - 8000 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	Matched Antibody Pair kit is a titrated unlabeled capture antibody, a titrated biotin-labeled detector and a calibrated protein standard. The Matched Antibody Pair Kit can be used to quantify native and recombinant human MMP1.

Both capture and detector antibodies are rabbit monoclonal antibodies.

Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.

Additional protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® ([ab215083](#)) which uses the same antibody pair.

**To receive an electronic copy of the Certificate of Analysis, please send an email with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.**

Buffer information:  
 The capture antibody is glycerol free.  
 The detector antibody contains glycerol.

<b>Tested applications</b>	<b>Suitable for:</b> ELISA
<b>Platform</b>	Reagents

## Properties

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

Components	10 x 96 tests	2 x 96 tests
Human MMP1 Capture Antibody	1 x 100µg	1 x 20µg
Human MMP1 Detector Antibody	1 x 25µg	1 x 5µg
Human MMP1 Lyophilized Protein	1 vial	1 vial

**Function** Cleaves collagens of types I, II, and III at one site in the helical domain. Also cleaves collagens of types VII and X. In case of HIV infection, interacts and cleaves the secreted viral Tat protein, leading to a decrease in neuronal Tat's mediated neurotoxicity.

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

**Domain** There are two distinct domains in this protein; the catalytic N-terminal, and the C-terminal which is involved in substrate specificity and in binding TIMP (tissue inhibitor of metalloproteinases). 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.

**Post-translational modifications** Undergoes autolytic cleavage to two major forms (22 kDa and 27 kDa). A minor form (25 kDa) is the glycosylated form of the 22 kDa form. The 27 kDa form has no activity while the 22/25 kDa form can act as activator for collagenase.

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

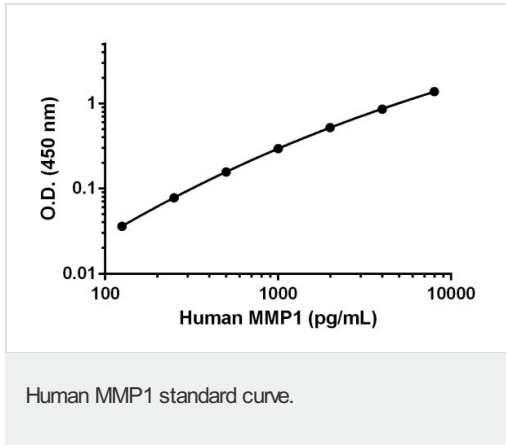
## Applications

Our [Abpromise guarantee](#) covers the use of **ab218184** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

## Images



Standard calibration curve. Background subtracted values are graphed.

**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