

# Human MMP9 Antibody Pair - BSA and Azide free ab241787

Recombinant **RabMAb**

[1 Image](#)

### Overview

<b>Product name</b>	Human MMP9 Antibody Pair - BSA and Azide free
<b>Assay type</b>	ELISA set
<b>Range</b>	15.6 pg/ml - 1000 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	<p>The Antibody Pair can be used to quantify Human MMP9. BSA and Azide free antibody pairs include unconjugated capture and detector antibodies suitable for sandwich ELISAs. The antibodies are provided at an approximate concentration of 1 mg/ml as measured by the protein A280 method. The recommended antibody orientation is based on internal optimization for ELISA-based assays. Antibody orientation is assay dependent and needs to be optimized for each assay type. Both capture and detector antibodies are rabbit monoclonal antibodies delivering consistent, specific, and sensitive results.</p> <p>For additional information on the performance of the antibody pair, see the equivalent Matched Antibody Pair Kit (<a href="#">ab213741</a>), which uses the same antibodies. Please note that the range provided for the pairs is only an estimation based on the performance of the related product using the same antibody pair. Performance of the antibody pair will depend on the specific characteristics of your assay. We guarantee the product works in sandwich ELISA, but we do not guarantee the sensitivity or dynamic range of the antibody pair in your assay.</p>
<b>Tested applications</b>	<b>Suitable for:</b> Sandwich ELISA
<b>Platform</b>	Reagents

### Properties

<b>Storage instructions</b>	Store at +4°C. Please refer to protocols.
<b>Carrier free</b>	Yes

Components	10 x 96 tests
Human MMP9 Capture Antibody (unconjugated)	1 x 100µg

Components	10 x 96 tests
Human MMP9 Detector Antibody (unconjugated)	1 x 100µg

<b>Function</b>	May play an essential role in local proteolysis of the extracellular matrix and in leukocyte migration. Could play a role in bone osteoclastic resorption. Cleaves KiSS1 at a Gly-Leu bond. Cleaves type IV and type V collagen into large C-terminal three quarter fragments and shorter N-terminal one quarter fragments. Degrades fibronectin but not laminin or Pz-peptide.
<b>Tissue specificity</b>	Produced by normal alveolar macrophages and granulocytes.
<b>Involvement in disease</b>	Intervertebral disc disease Metaphyseal anadysplasia 2
<b>Sequence similarities</b>	Belongs to the peptidase M10A family. Contains 3 fibronectin type-II domains. Contains 4 hemopexin repeats.
<b>Domain</b>	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.
<b>Post-translational modifications</b>	Processing of the precursor yields different active forms of 64, 67 and 82 kDa. Sequentially processing by MMP3 yields the 82 kDa matrix metalloproteinase-9. N- and O-glycosylated.
<b>Cellular localization</b>	Secreted, extracellular space, extracellular matrix.

## Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab241787 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
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

## Images

Powered by  
recombinant antibodies



**Research with confidence**  
Consistent and reproducible results



**Long-term and scalable supply**  
Recombinant technology



**Success from the first experiment**  
Confirmed specificity



**Ethical standards compliant**  
Animal-free production

Sandwich ELISA - Human MMP9 Antibody Pair -  
BSA and Azide free (ab241787)

To learn more about the advantages of recombinant antibodies see [here](#).

**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