

# Human MMP9 Matched Antibody Pair Kit ab213741

重组

## 2 图像

### 概述

产品名称	人MMP9抗体对试剂盒
检测方法	Colorimetric
检测类型	ELISA set
灵敏度	3.77 pg/ml
范围	15.6 pg/ml - 1000 pg/ml
种属反应性	<b>与反应:</b> Human
产品概述	<p>Human MMP9 Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant human MMP9.</p> <p>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.</p> <p>Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our <a href="#">website</a>. An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs (<a href="#">ab210905</a>).</p>
平台	Microplate

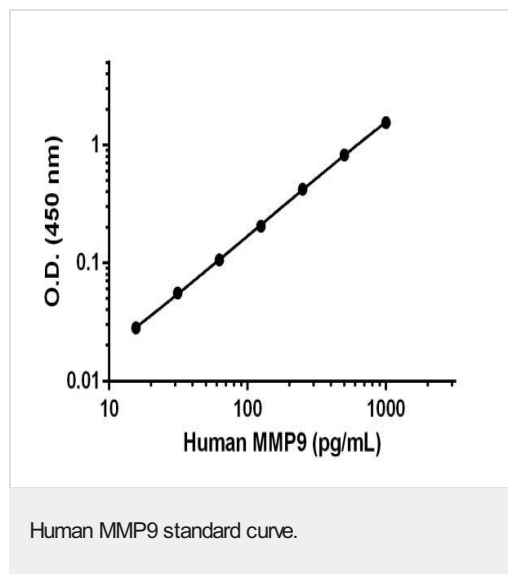
### 性能

**存放说明** Store at -20°C. Please refer to protocols.

组件	5 x 96 tests
Human MMP9 Capture Antibody	1 x 50µg
Human MMP9 Detector Antibody	1 x 12.5µg
Human MMP9 Lyophilized Protein	1 vial

<b>功能</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>组织特异性</b>	Produced by normal alveolar macrophages and granulocytes.
<b>疾病相关</b>	Intervertebral disc disease Metaphyseal anadysplasia 2
<b>序列相似性</b>	Belongs to the peptidase M10A family. Contains 3 fibronectin type-II domains. Contains 4 hemopexin repeats.
<b>结构域</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>翻译后修饰</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>细胞定位</b>	Secreted, extracellular space, extracellular matrix.

## 图片



Standard calibration curve. Background subtracted values are graphed. Data provided for demonstration purposes only.

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 Matched Antibody  
Pair Kit (ab213741)

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.cn/abpromise> or contact our technical team.

### Terms and conditions

---

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