

Mouse proMMP9 ELISA Kit ab100732

5 References 1 图像

概述		
产品名称	小鼠proMMP9 ELISA试剂盒	
检测方法	Colorimetric	
样品类型	Cell culture supernatant, Serum, Plasma	
检测类型	Sandwich (quantitative)	
灵敏度	< 50 pg/ml	
范围	74.07 pg/ml - 18000 pg/ml	
回收率	100 %	
		特定样本回收率
样品类型		平均 %
Cell culture supernatant		109.7
Serum		114.4
Plasma		112
		范围
		96% - 118%
		97% - 132%
		101% - 121%
实验步骤	Multiple steps standard assay	
种属反应性	与反应: Mouse	
产品概述	<p>proMMP9 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab100732) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Mouse proMMP9 in serum, plasma and cell culture supernatants.</p> <p>This assay employs an antibody specific for mouse proMMP9 coated on a 96-well plate. Standards and samples are pipetted into the wells and proMMP9 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse proMMP9 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 proMMP9 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>	
平台	Microplate	

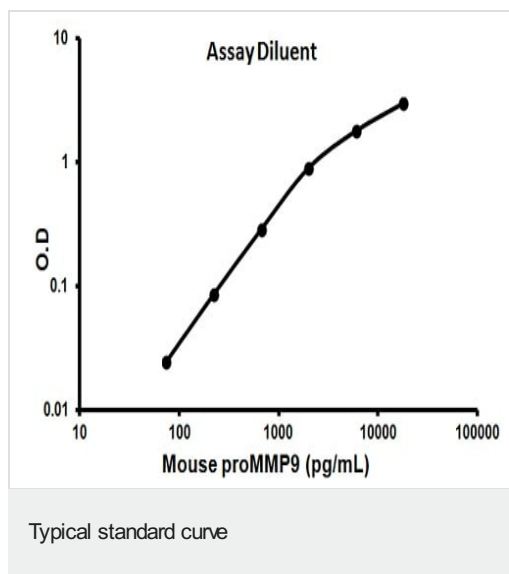
性能

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-Mouse proMMP9	2 vials
proMMP9 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse proMMP9 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

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

图片



Representative standard curve using ab100732

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