

Mouse MIP2 ELISA Kit (CXCL2) ab204517

重组 SimpleStep ELISA

★★★★☆ [1 Abreviews](#) [5 References](#) [8 图像](#)

概述

产品名称 小鼠MIP2 ELISA试剂盒(CXCL2)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
overall	8			1.64%

批次间

样品	n	Mean	SD	CV%
overall	3			4.87%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 0.87 pg/ml

范围 3.13 pg/ml - 200 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Serum	108	103% - 113%
Cell culture media	102	98% - 107%
Hep Plasma	104	101% - 109%
EDTA Plasma	115	113% - 116%
Cit plasma	91	87% - 97%

检测时间 1h 30m

实验步骤 One step assay

种属反应性

与反应: Mouse

产品概述

Mouse MIP2 ELISA Kit (CXCL2) (ab204517) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of MIP2 (CXCL2) protein in cell culture supernatant, plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse MIP2 (CXCL2) with 0.87 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

Mouse Macrophage Inflammatory Protein-2 (MIP2), also known as C-X-C motif chemokine 2 (CXCL2), is a small cytokine belonging to the CXC chemokine family. MIP2 was originally identified as a heparin-binding protein, and has been shown to exhibit potent neutrophil chemotactic activity.

Mouse MIP2 c-DNA encodes a 100 amino acid residue precursor protein. The amino-terminal 27 amino acid residues are cleaved from this precursor to generate the mature mouse MIP2. Mouse MIP2 is 63% identical to Mouse KC (another mouse alpha chemokine), and mouse MIP2 is 60% identical to human GROβ and GROγ. Based on these protein sequence similarities, it is likely that mouse MIP2 and KC are homologs of the human GROα, β, and γ chemokines. However, since chemokines with protein sequence homology to human IL-8 have not been identified in mice, it has been suggested that the mouse MIP2 and KC are functional homologs of human IL-8 in mice. A putative mouse homolog of the human IL-8 receptor beta (IL-8 Rβ) has also been cloned. This receptor shows 71% identity to human IL-8 Rβ and 68% identity to human IL-8 Rα. Both mouse MIP2 and KC bind mouse IL-8 Rβ with high affinity.

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Mouse MIP2 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml

组件	1 x 96 tests
Mouse MIP2 Capture Antibody (Lyophilized)	1 vial
Mouse MIP2 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent 50BP	1 x 20ml
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

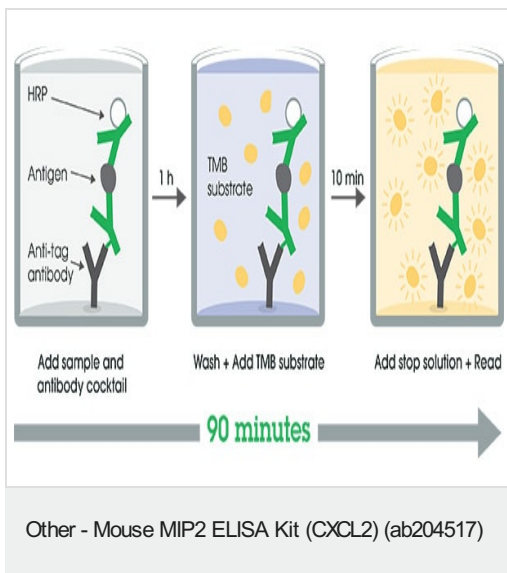
功能 Produced by activated monocytes and neutrophils and expressed at sites of inflammation. Hematopoietic chemokine, which, in vitro, suppresses hematopoietic progenitor cell proliferation. GRO-beta(5-73) shows a highly enhanced hematopoietic activity.

序列相似性 Belongs to the intercrine alpha (chemokine CxC) family.

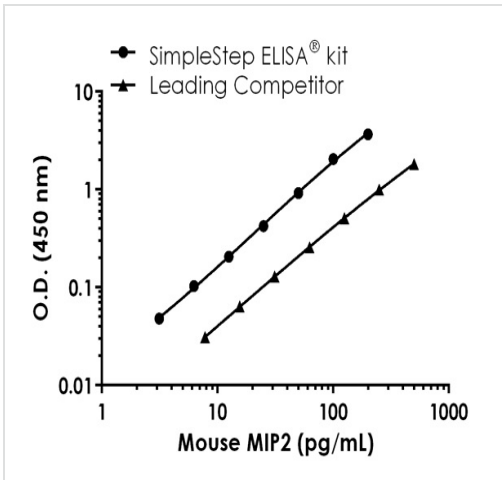
翻译后修饰 The N-terminal processed form GRO-beta(5-73) is produced by proteolytic cleavage after secretion from bone marrow stromal cells.

细胞定位 Secreted.

图片

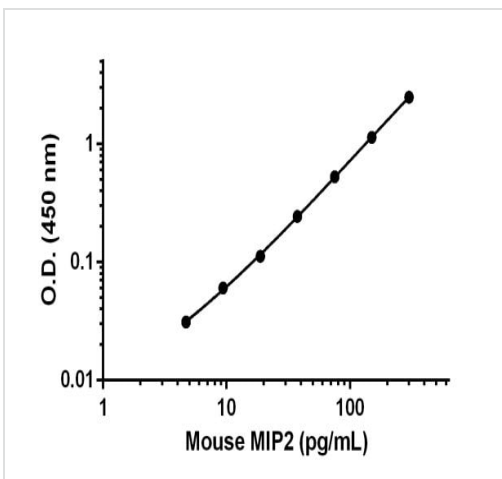


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



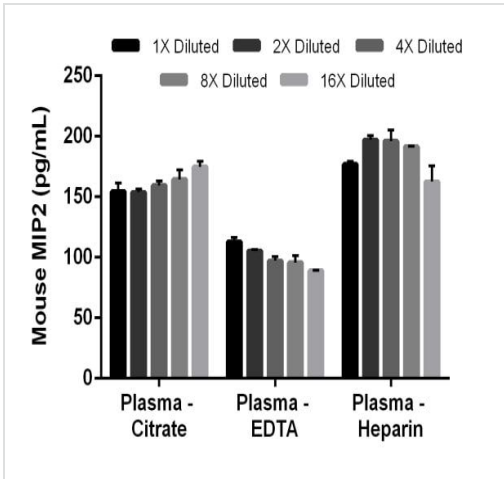
Mouse MIP2 standard curve comparison data.

Standard curve comparison between mouse MIP2 SimpleStep ELISA[®] kit and traditional ELISA kit from leading competitor. SimpleStep ELISA kit show comparable sensitivity.



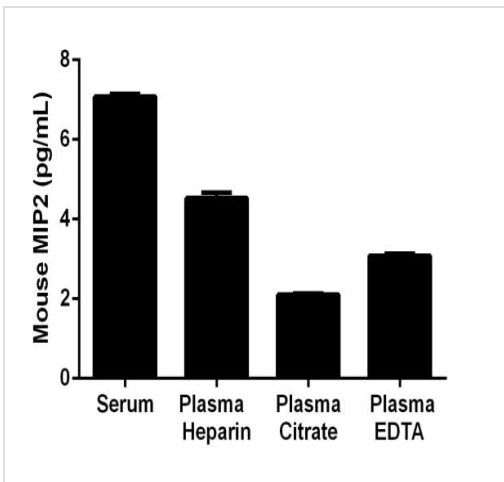
Linearity of dilution of MIP2 in serum and cell culture media.

Recombinant mouse MIP2 was spiked into 100% serum and diluted in a 2-fold dilution series in Sample Diluent 50BP. Recombinant mouse MIP2 was spiked into 10% cell culture media and diluted in a 2-fold dilution series in Sample Diluent NS. The concentrations of mouse MIP2 were measured in duplicate and interpolated from the mouse MIP2 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are graphed (mean +/- SD).



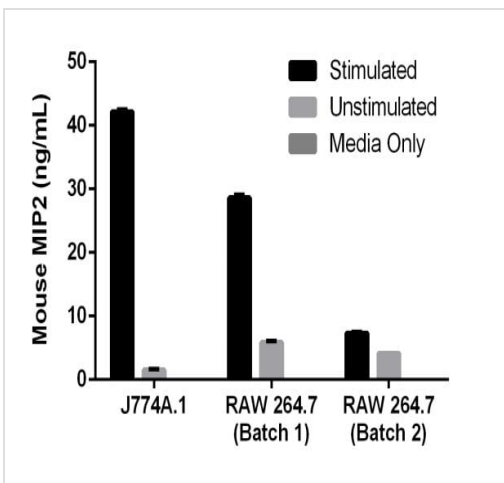
Linearity of dilution of spiked recombinant MIP2 in plasma.

Recombinant mouse MIP2 was spiked into 100% citrate plasma and 50% EDTA plasma and diluted in a 2-fold dilution series in Sample Diluent 25BP. Recombinant mouse MIP2 was spiked into 100% heparin plasma and diluted in a 2-fold dilution series in Sample Diluent 50BP. The concentrations of mouse MIP2 were measured in duplicate and interpolated from the mouse MIP2 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are graphed (mean +/- SD).



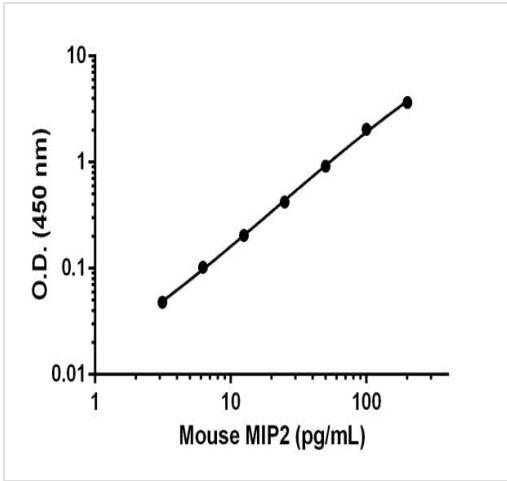
Mouse MIP2 expression is shown for serum and plasma samples.

Native MIP2 was measured in duplicate in 100% serum and 100% heparin plasma and concentrations interpolated from a standard curve diluted in Sample Diluent 50BP. Native MIP2 was measured in duplicate in 100% Citrate plasma and 50% EDTA plasma and concentrations interpolated from a standard curve diluted in Sample Diluent 25BP. The interpolated dilution factor corrected values are graphed (mean +/- SD).



Mouse MIP2 expression is shown for cultured media from two mouse cell lines.





J774A.1 cells were cultured in HGDMEM with 10% fetal calf serum, and 100 µg/mL of Kanamycin. During the exponential growth phase, J774A.1 cells were treated for 72 hours in the presence and absence of 1.5% PHA and 10 ng/mL of PMA. RAW 264.7 cells were cultured in HGDMEM with 10% fetal calf serum, 2 mM L-glutamine and 100 µg/mL Kanamycin. During the exponential growth phase, RAW264.7 cells were starved for 24 hours and treated in the presence and absence of 5 µg/mL of LPS (Batch #1) or 1% PHA (Batch #2). The concentrations of mouse MIP2 were interpolated from a standard curve diluted in Sample Diluent NS and corrected for sample dilution. The interpolated dilution factor corrected values are graphed (mean +/- SD).



Background-subtracted data values (mean +/- SD) are graphed.

Example of the mouse MIP2 standard curve in Sample Diluent NS.

Powered by recombinant antibodies

 <p>Research with confidence Consistent and reproducible results</p>	 <p>Long-term and scalable supply Recombinant technology</p>
 <p>Success from the first experiment Confirmed specificity</p>	 <p>Ethical standards compliant Animal-free production</p>

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

Sandwich ELISA - Mouse MIP2 ELISA Kit (CXCL2)
(ab204517)

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