

Rat MIP-1 beta ELISA Kit (CCL4) ab171336

1 References [画像数 1](#)

製品の概要

製品名 Rat MIP-1 beta ELISA Kit (CCL4)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
1	16	31pg/ml	1.42	= 4.6%
2	16	144pg/ml	9.21	= 6.4%
3	16	453pg/ml	24.31	= 5.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
1	24	31pg/ml	1.67	= 5.4%
2	24	148pg/ml	12.13	= 8.2%
3	24	419pg/ml	24.3	= 5.8%

サンプルの種類 Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 < 10 pg/ml

検出範囲 15.6 pg/ml - 1000 pg/ml

ステップ Multiple steps standard assay

種交差性 交差種: Rat

製品の概要 Abcam's rat MIP-1 beta (CCL4) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of rat MIP-1 beta in cell culture supernatants, serum and plasma (heparin, EDTA, citrate)

A MIP-1 beta specific rat monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection

polyclonal antibody from goat, specific for MIP-1 beta is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the rat MIP-1 beta amount of sample captured in plate.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	ラベル	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-rat MIP-1 beta Antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-rat MIP-1 beta antibody		1 x 100µl
Lyophilized recombinant rat MIP-1 beta standard		2 x 10ng
Plate Seal		4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

機能

Monokine with inflammatory and chemokinetic properties. Binds to CCR5. One of the major HIV-suppressive factors produced by CD8+ T-cells. Recombinant MIP-1-beta induces a dose-dependent inhibition of different strains of HIV-1, HIV-2, and simian immunodeficiency virus (SIV). The processed form MIP-1-beta(3-69) retains the abilities to induce down-modulation of surface expression of the chemokine receptor CCR5 and to inhibit the CCR5-mediated entry of HIV-1 in T-cells. MIP-1-beta(3-69) is also a ligand for CCR1 and CCR2 isoform B.

配列類似性

Belongs to the intercrine beta (chemokine CC) family.

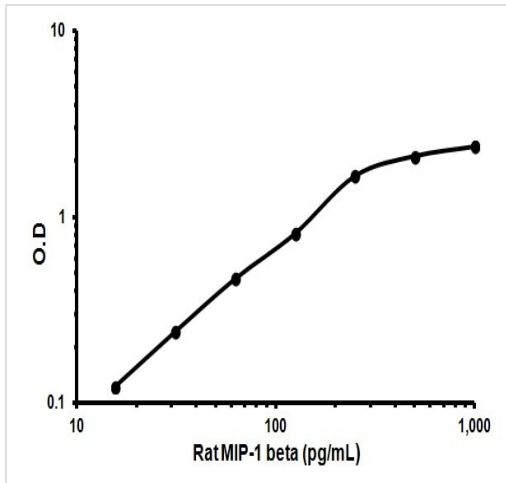
翻訳後修飾

N-terminal processed form MIP-1-beta(3-69) is produced by proteolytic cleavage after secretion from peripheral blood lymphocytes.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab171336.

Typical standard curve

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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