

Product datasheet

Rat MIP-1 beta ELISA Kit (CCL4) ab171336

[1 References](#) [1 Image](#)

Overview

Product name Rat MIP-1 beta ELISA Kit (CCL4)

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 10% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 10% |

Sample type Cell culture supernatant, Plasma, Platelets

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 15.6 pg/ml - 1000 pg/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat

Product overview

Abcam's rat MIP-1 beta (CCL4) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of mouse 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.

Tested applications **Suitable for:** Sandwich ELISA

Platform Microplate

Properties

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

| Components | Identifier | 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 130µl |
| Biotinylated anti-rat MIP-1 beta antibody | | 1 x 130µl |
| 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 |

Function 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.

Sequence similarities Belongs to the intercrine beta (chemokine CC) family.

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

Cellular localization Secreted.

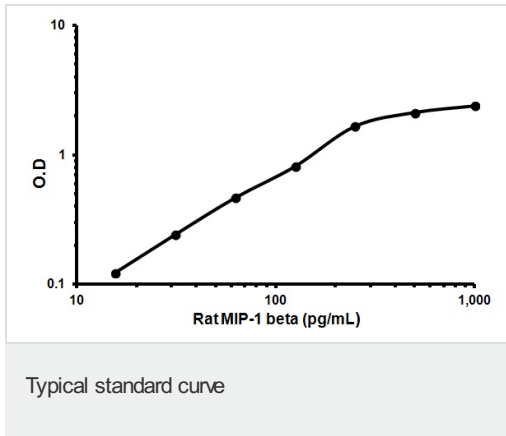
Applications

Our [Abpromise guarantee](#) covers the use of **ab171336** 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 |
|----------------|-----------|--|
| Sandwich ELISA | | Use at an assay dependent concentration. |

Images



Representative Standard Curve using ab171336.

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