

# Human MIP1 beta ELISA Kit (CCL4) ab100597

[2 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Human MIP1 beta ELISA Kit (CCL4)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 2.5 pg/ml
<b>Range</b>	4.1 pg/ml - 1000 pg/ml
<b>Recovery</b>	95 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	97.42	85% - 104%
Serum	95.64	82% - 102%
Plasma	94.32	83% - 103%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's MIP1 beta Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human MIP1 beta in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human MIP1 beta coated on a 96-well plate. Standards and samples are pipetted into the wells and MIP1 beta present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated antiHuman MIP-1 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 MIP1 beta bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Notes** Optimization may be required with urine samples.

This product may also recognise CC4L [C-C motif chemokine 4-like], a family member which

shares high homology with CCL4.

**Platform** Microplate

## Properties

---

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human MIP1 beta	2 vials
MIP1 beta Microplate (12 x 8 wells)	1 unit
Recombinant Human MIP1 beta Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**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.

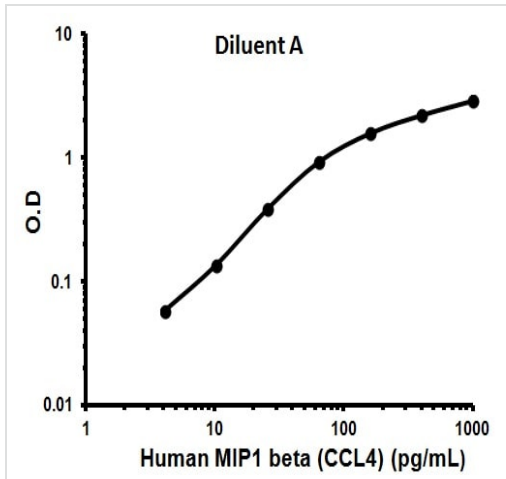
**Post-translational modifications** N-terminal processed form MIP-1-beta(3-69) is produced by proteolytic cleavage after secretion from peripheral blood lymphocytes.

**Cellular localization** Secreted.

---

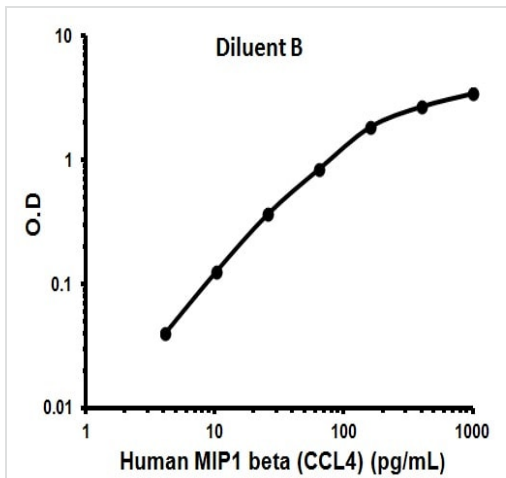
## Images

---



Representative Standard Curve using ab100597.

Typical Standard Curve



Representative Standard Curve using ab100597.

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.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