

Mouse MIP1α ELISA Kit (CCL3) αb100726

★☆☆☆☆ 1 Abreviews 1 References 1 Image

Overview

Product name	Mouse MIP1a ELISA Kit (CCL3)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 8 pg/ml
Range	4.1 pg/ml - 1000 pg/ml
Recovery	87 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	94.58	86% - 104%
Serum	88.58	83% - 99%
Plasma	87.94	82% - 98%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Mouse

Product overview Abcam’s MIP-1a Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse MIP-1a in serum, plasma, and cell culture supernatants.

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

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
240X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-Mouse MIP-1a	2 vials
MIP-1a Microplate (12 x 8 wells)	1 unit
Recombinant Mouse MIP-1a 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 CCR1, CCR4 and CCR5. One of the major HIV-suppressive factors produced by CD8+ T-cells. Recombinant MIP-1-alpha induces a dose-dependent inhibition of different strains of HIV-1, HIV-2, and simian immunodeficiency virus (SV).

Sequence similarities

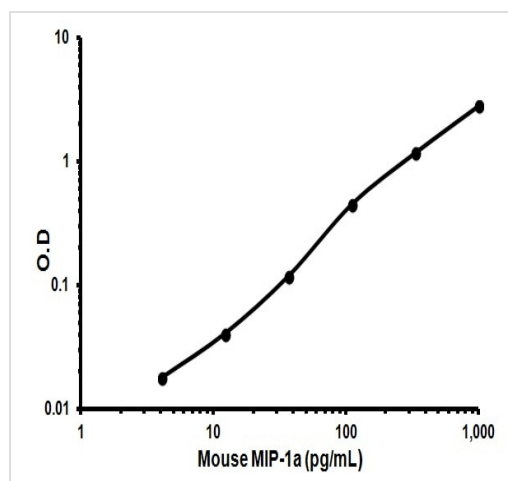
Belongs to the intercrine beta (chemokine CC) family.

Post-translational modifications

N-terminal processed form LD78-alpha(4-69) is produced by proteolytic cleavage after secretion from HTLV1-transformed T-cells.

Cellular localization

Secreted.

Images

Typical standard curve

Representative standard curve using ab100726

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