

Mouse MIP3 beta ELISA Kit (CCL19) ab100729

2 References 画像数 1

製品の概要

製品名	Mouse MIP3 beta ELISA Kit (CCL19)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 5 pg/ml
検出範囲	4.1 pg/ml - 1000 pg/ml
添加回収試験	100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	85.6	78% - 101%
Serum	107.7	82% - 132%
Plasma	119.5	83% - 140%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要

Abcam's MIP3 beta Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse MIP-3- beta in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse MIP3 beta coated on a 96- well plate. Standards and samples are pipetted into the wells and MIP3 beta present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse MIP3 beta 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 MIP3 beta bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

試験プラットフォーム Microplate

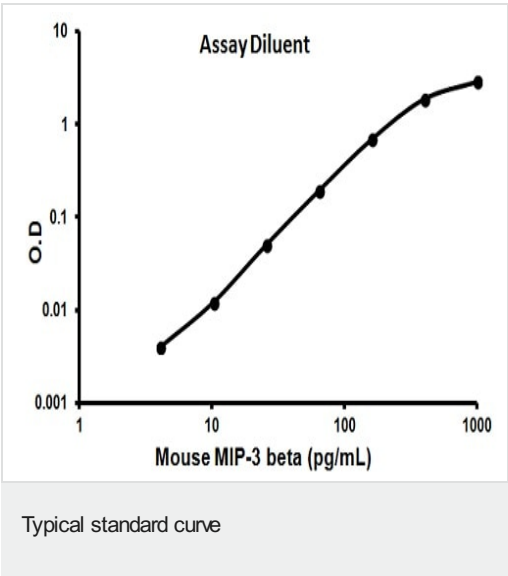
製品の特性

保存方法Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
260X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-Mouse MIP3 beta	2 vials
MIP3 beta Microplate (12 x 8 wells)	1 unit
Recombinant Mouse MIP3 beta Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能	May play a role not only in inflammatory and immunological responses but also in normal lymphocyte recirculation and homing. May play an important role in trafficking of T-cells in thymus, and T-cell and B-cell migration to secondary lymphoid organs. Specifically binds to chemokine receptor CCR7. Recombinant CCL19 shows potent chemotactic activity for T-cells and B-cells but not for granulocytes and monocytes.
組織特異性	Expressed at high levels in the lymph nodes, thymus and appendix. Intermediate levels seen in colon and trachea, while low levels found in spleen, small intestine, lung, kidney and stomach.
配列類似性	Belongs to the intercrine beta (chemokine CC) family.
細胞内局在	Secreted.

画像



Representative standard curve using ab100729

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