

Human MIP3b ELISA Kit (CCL19) ab100601

1 References 画像数 1

製品の概要

| | |
|---------|---|
| 製品名 | Human MIP3b 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 | < 2 pg/ml |
| 検出範囲 | 2.05 pg/ml - 500 pg/ml |
| 添加回収試験 | 78 % |

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 | 71.8 | 65% - 89% |
| Serum | 84.43 | 75% - 94% |
| Plasma | 79.88 | 71% - 90% |

| | |
|------------|--|
| ステップ | Multiple steps standard assay |
| 種交差性 | 交差種: Human |
| 製品の概要 | <p>Abcam's MIP3b Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human MIP3b in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human MIP3b coated on a 96- well plate. Standards and samples are pipetted into the wells and MIP3b present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human MIP3b 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 MIP3b bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p> |
| 特記事項 | Optimization may be required with urine samples. |
| 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 |
| 400X HRP-Streptavidin Concentrate | 1 x 200µl |
| 5X Assay Diluent | 1 x 15ml |
| Biotinylated anti-Human MIP3b | 2 vials |
| MIP3b Microplate (12 x 8 wells) | 1 unit |
| Recombinant Human MIP3b 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.

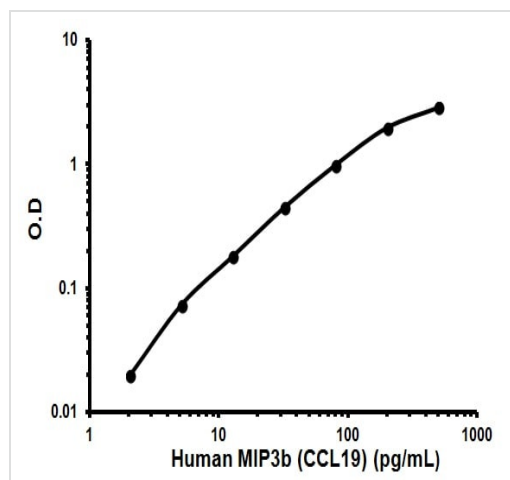
配列類似性

Belongs to the intercrine beta (chemokine CC) family.

細胞内局在

Secreted.

画像



Typical Standard Curve

Representative Standard Curve using ab100601.

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