

Human CCL23 ELISA Kit ab100611

2 References 画像数 2

製品の概要

製品名	Human CCL23 ELISA Kit
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	< 7 pg/ml
検出範囲	8.23 pg/ml - 6000 pg/ml
添加回収試験	93 %

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	93.48	81% - 105%
Serum	92.78	82% - 105%
Plasma	94.16	83% - 106%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's CCL23 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CCL23 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human CCL23 coated on a 96-well plate. Standards and samples are pipetted into the wells and CCL23 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CCL23 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 CCL23 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get higher sensitivity in only 90 minutes with Human CCL23 ELISA Kit ([ab216169](#)) from our SimpleStep ELISA® range.

特記事項 Optimisation 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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CCL23	2 vials
CCL23 Microplate (12 x 8 wells)	1 unit
Recombinant Human CCL23 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

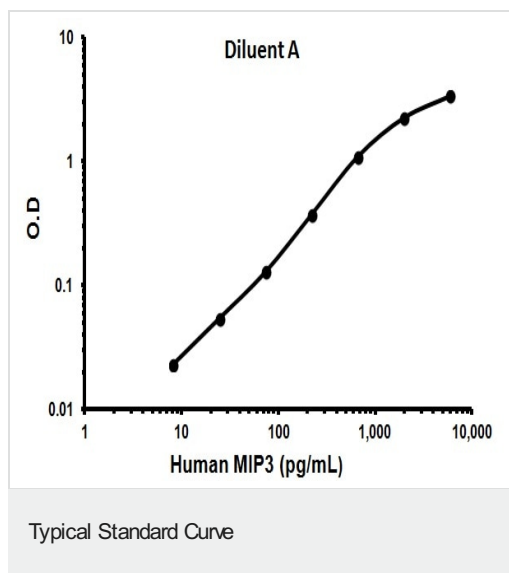
機能 Shows chemotactic activity for monocytes, resting T-lymphocytes, and neutrophils, but not for activated lymphocytes. Inhibits proliferation of myeloid progenitor cells in colony formation assays. This protein can bind heparin. Binds CCR1. CCL23(19-99), CCL23(22-99), CCL23(27-99), CCL23(30-99) are more potent chemoattractants than the small-inducible cytokine A23.

組織特異性 High levels in adult lung, liver, skeletal muscle and pancreas. Moderate levels in fetal liver, adult bone marrow and placenta. The short form is the major species and the longer form was detected only in very low abundance. CCL23(19-99), CCL23(22-99), CCL23(27-99), CCL23(30-99) are found in high levels in synovial fluids from rheumatoid patients.

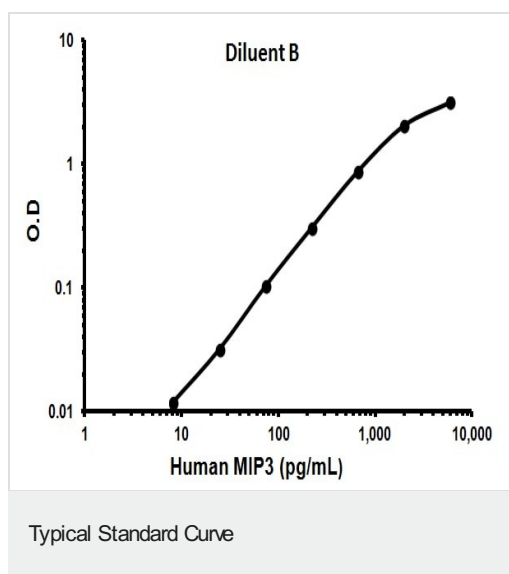
配列類似性 Belongs to the intercrine beta (chemokine CC) family.

細胞内局在 Secreted.

画像



Representative Standard Curve using ab100611.



Representative Standard Curve using ab100611.

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