

Human I-309 ELISA Kit (CCL1) ab100536

★★★★★ [1 Abreviews](#) [1 References](#) [画像数 1](#)

製品の概要

製品名	Human I-309 ELISA Kit (CCL1)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 4 pg/ml
検出範囲	4.12 pg/ml - 3000 pg/ml
添加回収試験	95 %

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	79.9	70% - 103%
Plasma	97.65	88% - 110%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam’s I-309 (CCL-1) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human I-309 (CCL-1) in plasma and cell culture supernatants. (Human I-309 concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for Human I-309 coated on a 96-well plate. Standards and samples are pipetted into the wells and I-309 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human I-309 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 I-309 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect the endogenous Human I-309 in normal serum with ab100536, only in serum spiked with Human I-309.

特記事項 Optimization may be required with urine samples

## 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
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-Human I-309	2 vials
I-309 Microplate (12 x 8 wells)	1 unit
Recombinant Human I-309 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## 機能

Cytokine that is chemotactic for monocytes but not for neutrophils. Binds to CCR8.

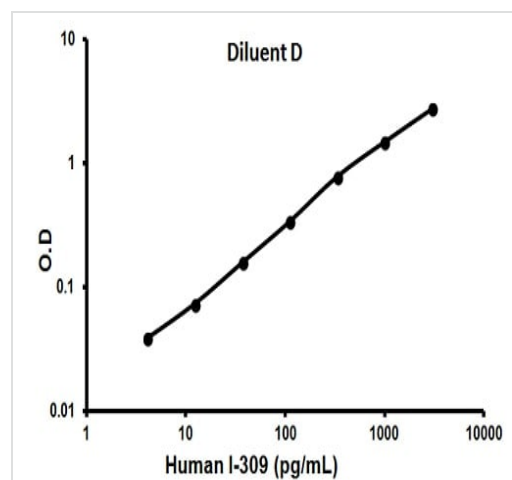
## 配列類似性

Belongs to the intercrine beta (chemokine CC) family.

## 細胞内局在

Secreted.

## 画像



Representative standard curve using ab100536

Typical standard curve

### **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