

### Human I-309 ELISA Kit (CCL1) ab100536

★★★★★ [1 Abreviews](#) [1 References](#) [1 图像](#)

#### 概述

产品名称	人I-309 ELISA试剂盒(CCL1)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 4 pg/ml
范围	4.12 pg/ml - 3000 pg/ml
回收率	95 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	79.9	70% - 103%
Plasma	97.65	88% - 110%

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** 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.

**说明** Optimization may be required with urine samples

平台

Microplate

性能

存放说明

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

组件	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

功能

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

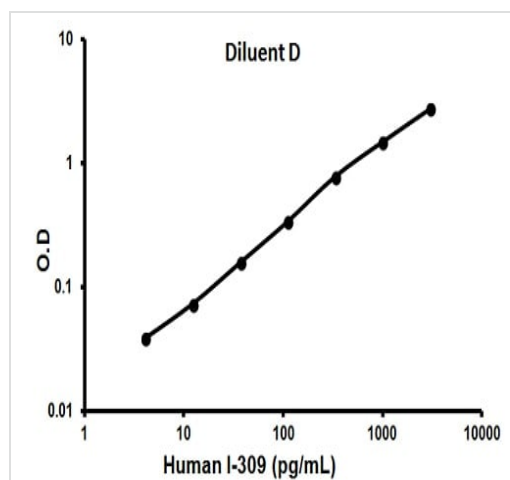
序列相似性

Belongs to the intercrine beta (chemokine CC) family.

细胞定位

Secreted.

图片



Representative standard curve using ab100536

Typical standard curve

**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.cn/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors