

Human CXCL9 ELISA Kit ab100595

★★★★★ [1 Abreviews](#) [2 References](#) [2 图像](#)

概述

产品名称	人CXCL9 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 20 pg/ml
范围	8.23 pg/ml - 6000 pg/ml
回收率	97 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	101.3	89% - 108%
Serum	94.56	82% - 102%
Plasma	96.32	84% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Abcam's CXCL9 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human CXCL9 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human CXCL9 coated on a 96-well plate. Standards and samples are pipetted into the wells and CXCL9 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human MIG 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 CXCL9 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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.

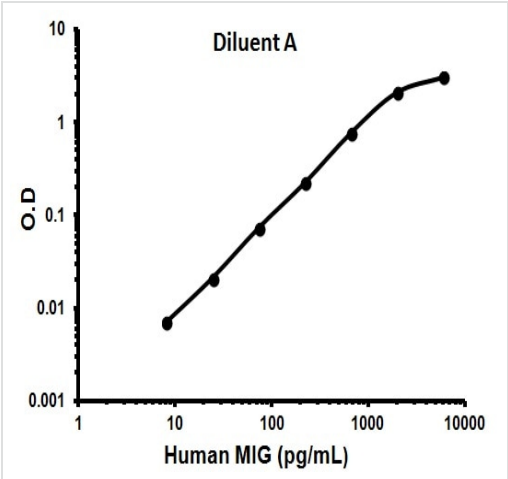
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CXCL9	2 vials
CXCL9 Microplate (12 x 8 wells)	1 unit
Recombinant Human CXCL9 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

功能Cytokine that affects the growth, movement, or activation state of cells that participate in immune and inflammatory response. Chemotactic for activated T-cells. Binds to CXCR3.

序列相似性Belongs to the intercrine alpha (chemokine CxC) family.

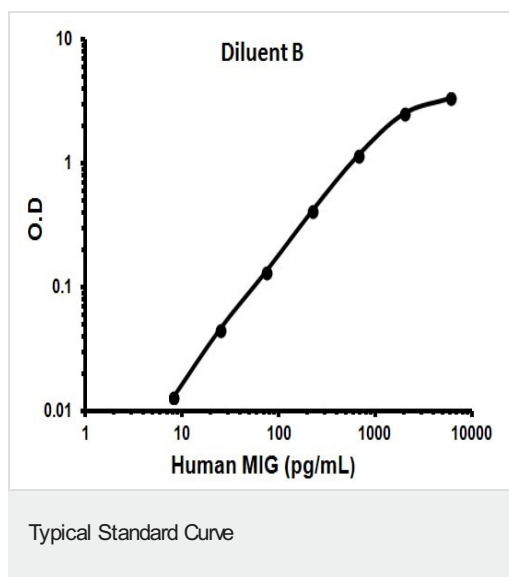
细胞定位Secreted.

图片



Typical Standard Curve

Representative Standard Curve using ab100595.



Representative Standard Curve using ab100595.

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