

Human Fractalkine ELISA Kit (CX3CL1) ab100522

2 References 2 图像

概述

| | |
|------|----------------------------------|
| 产品名称 | 人Fractalkine ELISA试剂盒(CX3CL1) |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Plasma |
| 检测类型 | Sandwich (quantitative) |
| 灵敏度 | < 300 pg/ml |
| 范围 | 0.25 ng/ml - 60 ng/ml |
| 回收率 | 93 % |

特定样本回收率

| 样品类型 | 平均% | 范围 |
|--------------------------|-------|------------|
| Cell culture supernatant | 94.47 | 83% - 102% |
| Plasma | 93.54 | 83% - 103% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's CX3CL1 (Fraktaline) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CX3CL1 in plasma and cell culture supernatants. (Human CX3CL1 concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for human CX3CL1 coated on a 96-well plate. Standards and samples are pipetted into the wells and CX3CL1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human CX3CL1 antibody is added. After washing away unbound biotinylated antibody, HRPconjugated 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 CX3CL1 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 endogenous Human Fractalkine in normal serum with ab100522, only in serum spiked with Human Fractalkine.

Get higher sensitivity in only 90 minutes with Human Fractalkine ELISA Kit (CX3CL1) ([ab192145](#)) from our SimpleStep ELISA[®] range.

说明 Optimization may be required with urine samples

平台 Microplate

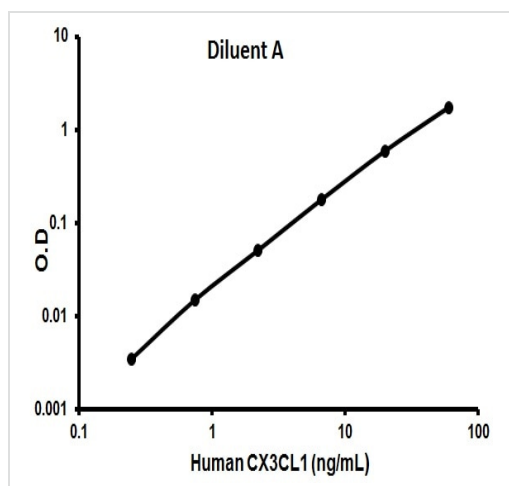
性能

存放说明 Store at -20°C. Please refer to protocols.

| 组件 | 1 x 96 tests |
|---|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 320X HRP-Streptavidin Concentrate | 1 x 200µl |
| 5X Assay Diluent B | 1 x 15ml |
| Assay Diluent A | 1 x 30ml |
| Biotinylated anti-Human CX3CL1 | 2 vials |
| CX3CL1 Microplate (12 x 8 well strips) | 1 unit |
| Recombinant Human CX3CL1 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

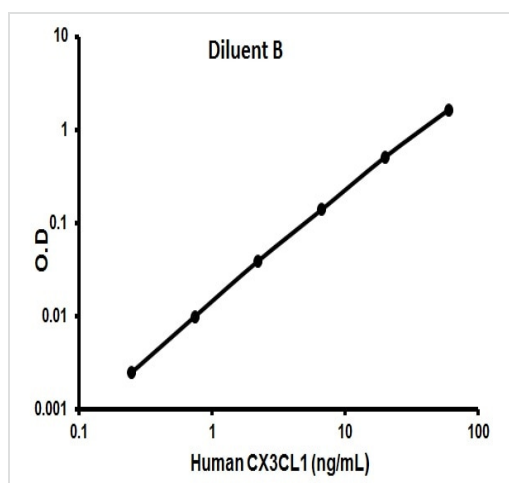
相关性 CX3CL1 is a member of the delta chemokine subfamily that contains a unique CX3C cysteine motif near the N-terminal. Unlike other known chemokines, it is a type 1 membrane protein containing a chemokine domain tethered on a long mucin-like stalk. CX3CL1, a leukocyte chemoattractant, is expressed in various tissues including the brain, heart, lung, kidney, skeletal muscle, and testis. The expression is reported to be up-regulated in endothelial cells and microglia by inflammatory signals. CX3CR1, a specific receptor for fractalkine, mediates both leukocyte migration and adhesion.

图片



Typical Standard Curve

Representative Standard Curve using ab100522.



Typical Standard Curve

Representative Standard Curve using ab100522.

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