

Human PF4 ELISA Kit (CXCL4) ab100628

4 References 2 图像

概述

产品名称	人PF4 ELISA试剂盒(CXCL4)
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	20.58 pg/ml - 15000 pg/ml
回收率	100 %

特定样本回收率

样品类型	平均 %	范围
Cell culture supernatant	108.2	101% - 116%
Serum	122.4	110% - 130%
Plasma	119.6	111% - 127%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's PF4 (CXCL4) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human PF4 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human PF4 coated on a 96-well plate. Standards and samples are pipetted into the wells and PF4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human PF4 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 PF4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human PF4 ELISA Kit (CXCL4) ([ab189573](#)) 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
500X HRP-Streptavidin	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Human PF4	2 units
PF4 Microplate (12 x 8 wells)	1 unit
Recombinant Human PF4 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

功能

Released during platelet aggregation. Neutralizes the anticoagulant effect of heparin because it binds more strongly to heparin than to the chondroitin-4-sulfate chains of the carrier molecule. Chemotactic for neutrophils and monocytes. Inhibits endothelial cell proliferation, the short form is a more potent inhibitor than the longer form.

序列相似性

Belongs to the intercrine alpha (chemokine CxC) family.

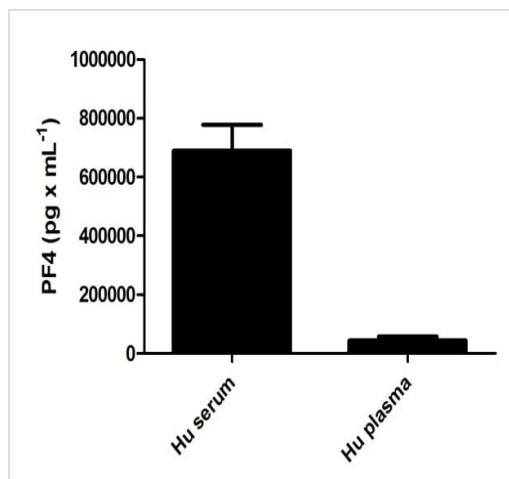
翻译后修饰

Binds non-covalently to a proteoglycan molecule.

细胞定位

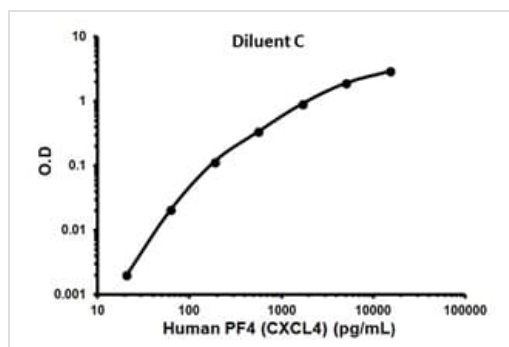
Secreted.

图片



Human PF4 measured in biological fluids showing quantity (pg) per mL of tested sample. Samples were diluted 200-800 fold.

Sandwich ELISA - PF4 (CXCL4) Human ELISA Kit  
(ab100628)



Representative Standard Curve using ab100628.

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