

Human PF4 ELISA Kit (CXCL4) ab100628

4 References 画像数 2

製品の概要

製品名	Human PF4 ELISA Kit (CXCL4)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 20 pg/ml
検出範囲	20.58 pg/ml - 15000 pg/ml
添加回収試験	100 %

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	108.2	101% - 116%
Serum	122.4	110% - 130%
Plasma	119.6	111% - 127%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

特記事項	Optimization may be required with urine samples.
試験プラットフォーム	Microplate

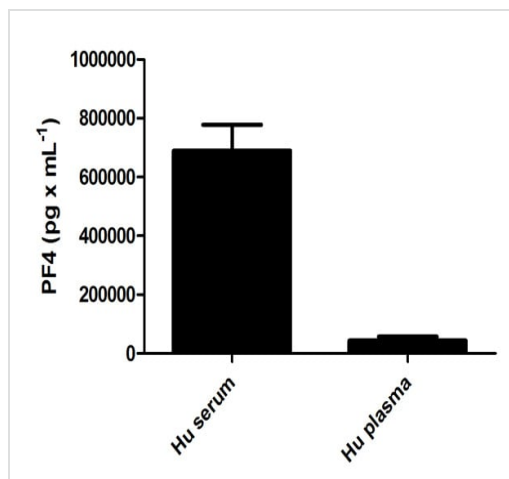
製品の特性

保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
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

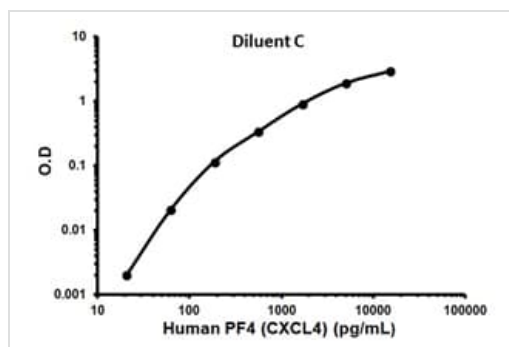
機能	Released during platelet aggregation. Neutralizes the anticoagulant effect of heparin because it binds more strongly to heparin than to the chondroitin-4-sulfate chains of the carrier molecule. Chemotactic for neutrophils and monocytes. Inhibits endothelial cell proliferation, the short form is a more potent inhibitor than the longer form.
配列類似性	Belongs to the intercrine alpha (chemokine CxC) family.
翻訳後修飾	Binds non-covalently to a proteoglycan molecule.
細胞内局在	Secreted.

画像



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.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