

Human PF4 ELISA Kit (CXCL4) ab189573

重组 SimpleStep ELISA

7 References 7 图像

概述

产品名称	人PF4 ELISA试剂盒(CXCL4)			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Overall	8		CV%
	批次间			
	样品	n	Mean	SD
	Overall	3		CV%
样品类型	Cell culture supernatant, Serum, Cell culture media, EDTA Plasma, Cit plasma			
检测类型	Sandwich (quantitative)			
灵敏度	24.29 pg/ml			
范围	78.13 pg/ml - 5000 pg/ml			
回收率	特定样本回收率			
	样品类型	平均%	范围	
	Cell culture supernatant	108	100% - 115%	
	Serum	93	89% - 99%	
	Cell culture media	87	85% - 90%	
	EDTA Plasma	103	97% - 115%	
	Cit plasma	92	83% - 100%	
检测时间	1h 30m			
实验步骤	One step assay			

种属反应性

与反应: Human

产品概述

As of November 23, 2020, Human PF4 (CXCL4) ELISA kit has been re-developed. We have identified new recombinant monoclonal antibodies to provide improved performance and consistency.

Human PF4 ELISA Kit (CXCL4) (ab189573) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of PF4 (CXCL4) protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human PF4 (CXCL4) with 24.29 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

Platelet factor-4 (PF4/CXCL4) is a 70-amino acid protein that is released from the alpha-granules of activated platelets whose major physiologic role appears to be inhibiting local antithrombin III activity and promoting coagulation. PF4 (CXCL4) has a role in inflammation and wound repair by acting as a strong chemoattractant for neutrophils, monocytes, and fibroblasts.

平台

Microplate

性能

存放说明

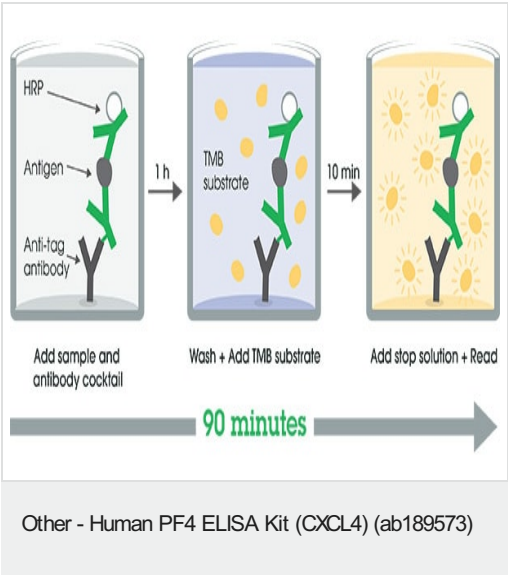
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests	1 x 96 tests
10X Human PF4 Capture Antibody	1 x 600µl	1 x 600µl
10X Human PF4 Detector Antibody	1 x 600µl	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 20ml
Antibody Diluent CPI2	1 x 6ml	1 x 6ml
Human PF4 Lyophilized Recombinant Protein (ab9562)	2 vials	2 vials
Plate Seals	1 unit	1 unit

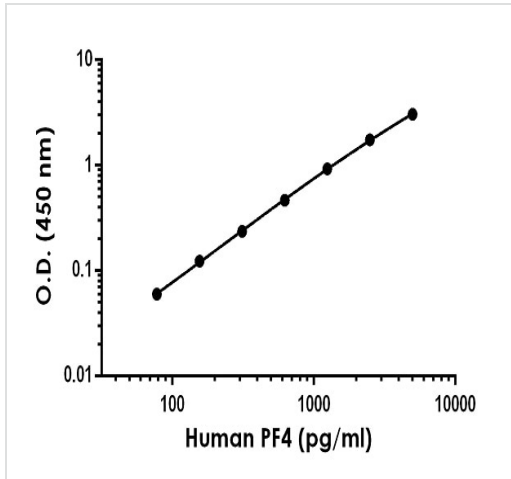
组件	1 x 96 tests	1 x 96 tests
Sample Diluent NS (ab193972)	1 x 50ml	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	1 unit
Stop Solution	1 x 12ml	1 x 12ml
TMB Development Solution	1 x 12ml	1 x 12ml

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

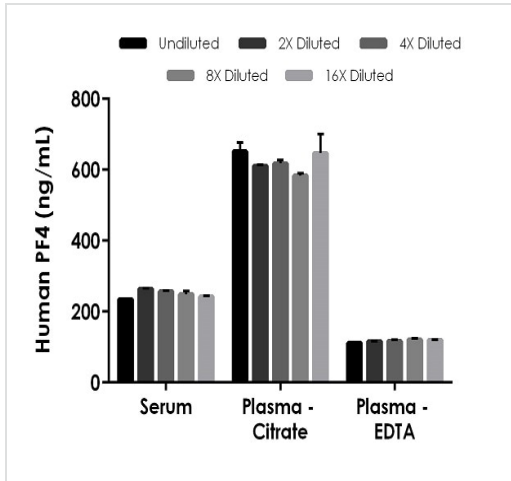


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



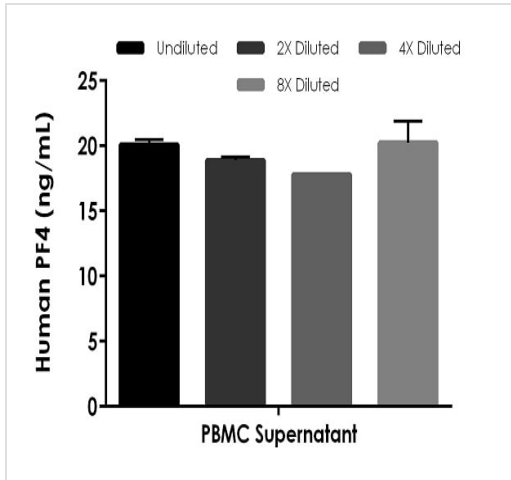
Raw data values are shown in the table. Background-subtracted data values (mean \pm SD) are graphed.

Example of human PF4 standard curve in Sample Diluent NS.



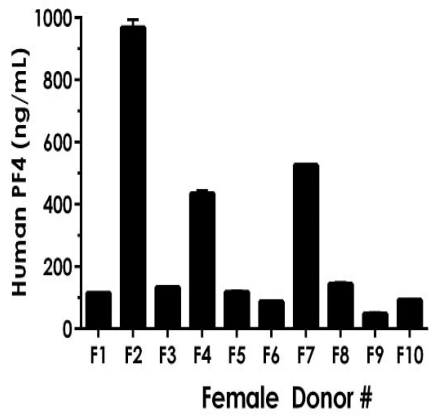
The concentrations of PF4 were measured in duplicates, interpolated from the PF4 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 2%, plasma (citrate) 0.5%, and plasma (EDTA) 4%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean PF4 concentration was determined to be 249.53 ng/mL in serum, 622.09 ng/mL in plasma (citrate), and 117.28 ng/mL in plasma (EDTA).

Interpolated concentrations of native PF4 in human serum, plasma and cell culture supernatant samples.



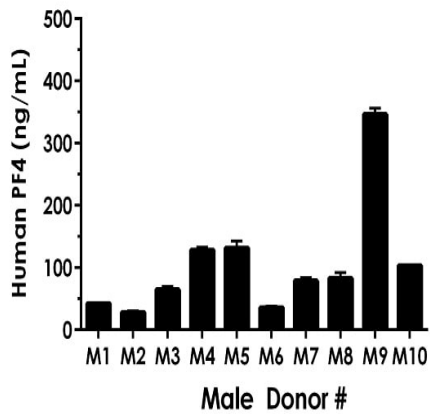
The concentrations of PF4 were measured in duplicates, interpolated from the PF4 standard curves and corrected for sample dilution. Undiluted samples are as follows: PBMC cell culture supernatant 12.5%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean PF4 concentration was determined to be 19.25 ng/mL in PBMC cell culture supernatant.

Interpolated concentrations of native PF4 in human cell culture supernatant samples.



Interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean PF4 concentration was determined to be 267.15 ng/mL with a range of 47.05 – 986.18 ng/mL.

0.5% Serum from ten individual healthy human female donors was measured in duplicate.



Interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean PF4 concentration was determined to be 104.38 ng/mL with a range of 27.73 – 353.63 ng/mL.

0.5% Serum from ten individual healthy human male donors was measured in duplicate.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human PF4 ELISA Kit (CXCL4)
(ab189573)

To learn more about the advantages of recombinant antibodies see [here](#).

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