

ab100628 – PF4 (CXCL4) Human ELISA Kit

Instructions for Use

For the quantitative measurement of Human PF4 (CXCL4) in serum, plasma, and cell culture supernatants.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

INTRODUCTION

1. BACKGROUND	2
2. ASSAY SUMMARY	3

GENERAL INFORMATION

3. PRECAUTIONS	4
4. STORAGE AND STABILITY	4
5. MATERIALS SUPPLIED	4
6. MATERIALS REQUIRED, NOT SUPPLIED	5
7. LIMITATIONS	5
8. TECHNICAL HINTS	6

ASSAY PREPARATION

9. REAGENT PREPARATION	7
10. STANDARD PREPARATIONS	8
11. SAMPLE PREPARATION	10
12. PLATE PREPARATION	10

ASSAY PROCEDURE

13. ASSAY PROCEDURE	11
---------------------	----

DATA ANALYSIS

14. CALCULATIONS	12
15. TYPICAL DATA	13
16. CALIBRATION	14
17. TYPICAL SAMPLE VALUES	14
18. ASSAY SPECIFICITY	15

RESOURCES

19. TROUBLESHOOTING	16
20. NOTES	17

1. BACKGROUND

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.

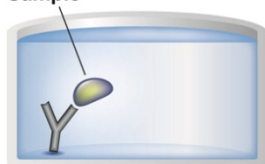
2. ASSAY SUMMARY

Primary Capture Antibody



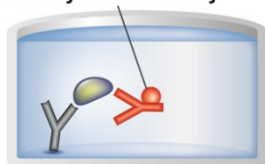
Prepare all reagents, samples and standards as instructed.

Sample



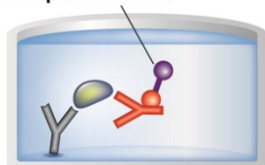
Add standard or sample to each well used. Incubate at room temperature

Biotinylated Antibody



Add prepared biotin antibody to each well. Incubate at room temperature.

Streptavidin-HRP



Add prepared Streptavidin solution. Incubate at room temperature.

Substrate Colored Product



Add TMB One-Step Development Solution to each well. Incubate at room temperature. Add Stop Solution to each well. Read at 450nm immediately.

3. PRECAUTIONS

Please read these instructions carefully prior to beginning the assay.

Modifications to the kit components or procedures may result in loss of performance.

4. STORAGE AND STABILITY

Store kit at -20°C immediately upon receipt.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in section 9. Reagent Preparation.

5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
PF4 Microplate (12 x 8 wells)	96 wells	-20°C
20X Wash Buffer Concentrate	25 mL	-20°C
Assay Diluent C	30 mL	-20°C
5X Assay Diluent B	15 mL	-20°C
Biotinylated anti-Human PF4	2 vials	-20°C
Recombinant Human PF4 Standard	2 vials	-20°C
500X HRP-Streptavidin Concentrate	200 µL	-20°C
TMB One-Step Substrate Reagent	12 mL	-20°C
Stop Solution	8mL	-20°C

6. MATERIALS REQUIRED, NOT SUPPLIED

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 2 μ L to 1 mL volumes.
- Adjustable 1-25 mL pipettes for reagent preparation.
- 100 mL and 1 liter graduated cylinders.
- Absorbent paper.
- Distilled or deionized water.
- Log-log graph paper or computer and software for ELISA data analysis.
- Tubes to prepare standard or sample dilutions.

7. LIMITATIONS

- Do not mix or substitute reagents or materials from other kit lots or vendors.

8. TECHNICAL HINTS

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps.
- When preparing your standards, it is very critical to briefly spin down the vial first. The powder may drop off from the cap when opening it if you do not spin down. Be sure to dissolve the powder thoroughly when reconstituting. After adding Assay Diluent to the vial, we recommend inverting the tube a few times, then flick the tube a few times, and then spin it down; repeat this procedure 3-4 times. This is a technique we find very effective for thoroughly mixing the standard without too much mechanical force.
- Do not vortex the standard during reconstitution, as this will destabilize the protein.
- Once your standard has been reconstituted, it should be used right away or else frozen for later use.
- Keep the standard dilutions on ice while during preparation, but the ELISA procedure should be done at room temperature.
- Be sure to discard the working standard dilutions after use – they do not store well.
- **This kit is sold based on number of tests. A ‘test’ simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.**

9. REAGENT PREPARATION

Equilibrate all reagents to room temperature (18-25°C) prior to use.

9.1 1X Assay Diluent B

5X Assay Diluent B should be diluted 5-fold with deionized or distilled water before use.

9.2 1X Wash Solution

If the 20X Wash Concentrate contains visible crystals, equilibrate to room temperature and mix gently until dissolved. Dilute 20 mL of 20X Wash Solution Concentrate into deionized or distilled water to yield 400 mL of 1X Wash Solution.

9.3 1X Biotinylated PF4 Detection Antibody

Briefly spin the Biotinylated anti-Human PF4 vial before use. Add 100 μ L of 1X Assay Diluent B into the vial to prepare a detection antibody concentrate. Pipette up and down to mix gently (the concentrate can either be stored at 4°C for 5 days or aliquoted and frozen at -20°C for 2 months). The detection antibody concentrate must be diluted 80-fold with 1X Assay Diluent B prior to use in the Assay Procedure.

9.4 1X HRP-Streptavidin Solution

Briefly spin the 500X HRP-Streptavidin concentrate vial before use. HRP-Streptavidin concentrate must be diluted 500-fold with 1X Assay Diluent B prior to use in the Assay Procedure.

For example: Briefly spin the vial and pipette up and down to mix gently. Add 20 μ L of 500X HRP-Streptavidin concentrate into a tube with 10 mL 1X Assay Diluent B to prepare a final 500-fold diluted 1X HRP-Streptavidin solution. Mix well.

10. STANDARD PREPARATIONS

- Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.
- Standard (recombinant protein) should be stored at -20°C or -80°C (recommended at -80°C) after reconstitution.

10.1 Briefly spin the vial of PF4 Standard. Prepare a 140 ng/mL PF4 **Stock Standard** by adding 400 µL Assay Diluent C.

10.2 Dissolve the powder thoroughly by gentle mixing.

10.3 Label tubes #1-7.

10.4 Prepare **Standard #1** by adding 75 µL 140 ng/mL **Stock Standard** to 625 µL Assay Diluent C into tube 1#. Mix thoroughly and gently.

10.5 Pipette 400 µL Assay Diluent C into each tube.

10.6 Prepare **Standard #2** by transferring 200 µL from tube #1 to #2, mix thoroughly.

10.7 Prepare **Standard #3** by transferring 200 µL from tube #2 to #3, mix thoroughly.

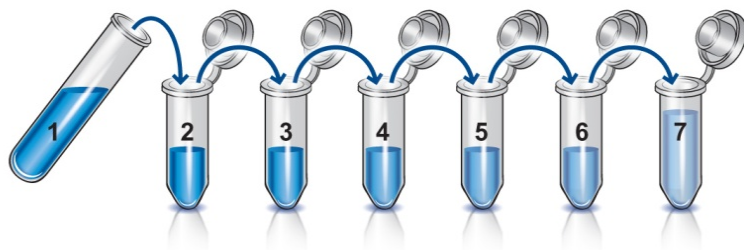
10.1 Using the table below as a guide, prepare further serial dilutions.

10.2 Assay Diluent C serves as the zero standard (0 pg/mL).

ASSAY PREPARATION

Standard Dilution Preparation Table

Standard #	Volume to Dilute (μL)	Diluent (μL)	Total Volume (μL)	Starting Conc. (pg/mL)	Final Conc. (pg/mL)
1	75	625	700	140,000	15,000
2	200	400	600	15,000	5,000
3	200	400	600	5,000	1,667
4	200	400	600	1,667	555.6
5	200	400	600	555.6	185.2
6	200	400	600	185.2	61.73
7	200	400	600	61.73	20.58
8	0	400	400	0	0



11. SAMPLE PREPARATION

General Sample Information:

- If your samples need to be diluted, Assay Diluent C should be used for dilution of serum, plasma and culture supernatants.
- Suggested dilution for normal serum/plasma: 200-2,000 fold.
- Please note that levels of the target protein may vary between different specimens. Optimal dilution factors for each sample must be determined by the investigator.

12. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused well strips should be returned to the plate packet and stored at 4°C.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

13. ASSAY PROCEDURE

- **Equilibrate all materials and prepared reagents to room temperature (18 - 25°C) prior to use.**
- **It is recommended to assay all standards, controls and samples in duplicate.**

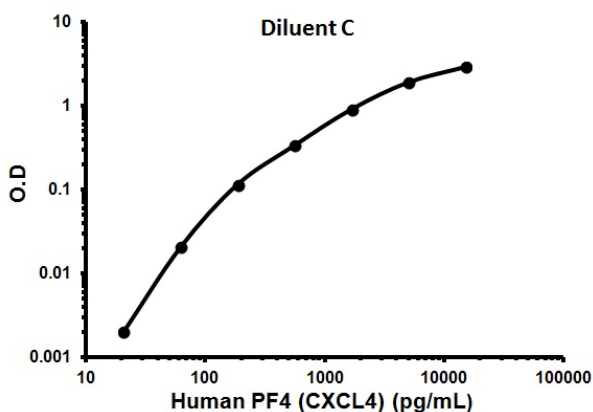
- 13.1 Add 100 µL of each standard (see Standard Preparation section 10) and sample into appropriate wells. Cover well and incubate for 2.5 hours at room temperature or over night at 4°C with gentle shaking.
- 13.2 Discard the solution and wash 4 times with 1X Wash Solution. Wash by filling each well with 1X Wash Solution (300 µL) using a multi-channel Pipette or autowasher. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining 1X Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
- 13.3 Add 100 µL of 1X Biotinylated PF4 Detection Antibody (Reagent Preparation section 9.3) to each well. Incubate for 1 hour at room temperature with gentle shaking.
- 13.4 Discard the solution. Repeat the wash as in step 13.2.
- 13.5 Add 100 µL of 1X HRP-Streptavidin solution (see Standard Preparation section 10) to each well. Incubate for 45 minutes at room temperature with gentle shaking.
- 13.6 Discard the solution. Repeat the wash as in step 13.2.
- 13.7 Add 100 µL of TMB One-Step Substrate Reagent to each well. Incubate for 30 minutes at room temperature in the dark with gentle shaking.
- 13.8 Add 50 µL of Stop Solution to each well. Read at 450 nm immediately.

14. CALCULATIONS

Calculate the mean absorbance for each set of duplicate standards, controls and samples, and subtract the average zero standard optical density. Plot the standard curve on log-log graph paper, with standard concentration on the x-axis and absorbance on the y-axis. Draw the best-fit straight line through the standard points.

15. TYPICAL DATA

TYPICAL STANDARD CURVE – Data provided for **demonstration purposes only**. A new standard curve must be generated for each assay performed.



Conc. (pg/mL)	O.D.
	Assay Diluent C
20.58	0.002
61.73	0.021
185.2	0.115
555.6	0.336
1,667	0.907
5,000	1.903
15,000	2.897

16. CALIBRATION

This immunoassay is calibrated against a highly purified Human PF4. The NIBSC/WHO unclassified purified Human PF4 preparation 83/505 was evaluated in this kit.

The dose response curve of the unclassified standard 83/505 parallels the PF4 standard curve. To convert sample values obtained with PF4 (CXCL4) Human ELISA Kit (ab100628) to approximate NIBSC Units/mL, use the equation below.

NIBSC (83/505) approximate value (Units/mL) = $2.57 \times \text{PF4 value (ng/mL)}$.

17. TYPICAL SAMPLE VALUES

SENSITIVITY –

The minimum detectable dose of PF4 is typically less than 20 pg/mL.

RECOVERY –

Recovery was determined by spiking various levels of Human PF4 into Human serum, plasma and cell culture media. Mean recoveries are as follows:

Sample Type	Average % Recovery	Range (%)
Serum	122.4	110-130
Plasma	119.6	111-127
Cell Culture Media	108.2	101-116

LINEARITY OF DILUTION -

Serum Dilution	Average % Expected Value	Range (%)
1:2	121.5	112-30
1:4	77.28	70-89

Plasma Dilution	Average % Expected Value	Range (%)
1:2	125.2	114-132
1:4	82.84	74-91

Cell Culture Media Dilution	Average % Expected Value	Range (%)
1:2	130.2	116-138
1:4	86.55	78-95

PRECISION –

	Intra-Assay	Inter-Assay
CV (%)	<10%	<12%

18. ASSAY SPECIFICITY

Cross Reactivity: This ELISA kit shows no cross-reactivity with any of the cytokines tested (e.g., Human Angiogenin, BDNF, BLC, ENA-78, FGF4, IL-1 α , IL-1 β , IL-2, IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-11, IL-12 p70, IL-12 p40, IL-13, IL-15, IL-309, IP-10, G-CSF, GM-CSF, IFN- γ , Leptin, MCP-1, MCP-3, MDC, MIP-1 α , MIP-1 β , MIP-1 δ , MMP-1,-2,-3,-10, PARC, RANTES, SCF, TARC, TGF- β , TIMP-1, TIMP-2, TNF- α , TNF- β , TPO, VEGF).

19. TROUBLESHOOTING

Problem	Cause	Solution
Poor standard curve	Inaccurate pipetting	Check pipettes
	Improper standards dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times; change to overnight standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	Store the reconstituted protein at -80°C, all other assay components 4°C. Keep substrate solution protected from light.
	Stop solution	Stop solution should be added to each well before measure.

20. NOTES

UK, EU and ROW

Email: technical@abcam.com | Tel: +44-(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com | Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel: 01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 400 921 0189 / +86 21 2070 0500

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp