

ab100781 – CXCL7 (PPBP) Rat ELISA Kit

Instructions for Use

For the quantitative measurement of Rat CXCL7 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. TYPICAL SAMPLE VALUES	14
17. ASSAY SPECIFICITY	15

RESOURCES

18. TROUBLESHOOTING	16
19. NOTES	17

1. BACKGROUND

Abcam's CXCL7 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat CXCL7 in serum, plasma, cell culture supernatants.

This assay employs an antibody specific for Rat CXCL7 coated on a 96-well plate. Standards and samples are pipetted into the wells and CXCL7 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat CXCL7 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 CXCL7 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

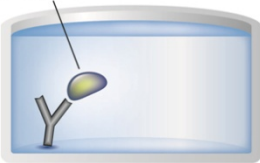
2. ASSAY SUMMARY

Primary Capture Antibody



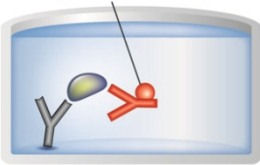
Prepare all reagents, samples and standards as instructed.

Sample



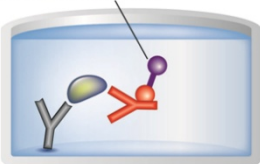
Add standard or sample to each well used. Incubate at room temperature

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)
CXCL7 Microplate (12 x 8 wells)	96 wells	-20°C
20X Wash Buffer Concentrate	25mL	-20°C
Recombinant Rat CXCL7 Standard	2 vials	-20°C
5X Assay Diluent	15 mL	-20°C
Biotinylated anti-Rat CXCL7	2 vials	-20°C
400X HRP-Streptavidin Concentrate	200 µL	-20°C
TMB One-Step Substrate Reagent	12 mL	-20°C
Stop Solution	8 mL	-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

5X Assay Diluent 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 Buffer Concentrate into deionized or distilled water to yield 400 mL of 1X Wash Buffer.

9.3 1X Biotinylated CXCL7 Detection Antibody

Briefly spin the Biotinylated- CXCL7 vial before use. Add 100 μ L of 1X Assay Diluent 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 prior to use in the Assay Procedure.

9.4 1X HRP-Streptavidin Solution

Briefly spin the 400X HRP-Streptavidin concentrate vial and pipette up and down to mix gently before use. HRP-Streptavidin concentrate must be diluted 400-fold with 1X Assay Diluent prior to use in the Assay Procedure.

For example: Briefly spin the vial and pipette up and down to mix gently. Add 30 μ L of 400X HRP-Streptavidin concentrate into a tube with 12 mL 1X Assay Diluent to prepare a final 400 fold diluted 1X HRP-Streptavidin solution (don't store the diluted solution for next day use). 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 CXCL7 Standard. Prepare the 100 ng/mL **Stock Standard** by adding 400 µL 1X Assay Diluent into the vial (see table below).

10.2 Ensure the powder is thoroughly dissolved by gentle mixing.

10.3 Label tubes #1-7.

10.4 Prepare **Standard #1** by adding 180 µL of the 100 ng/mL **Stock Standard**, to 180 µL of 1X Assay Diluent into tube #1. Mix thoroughly and gently.

10.5 Pipette 240µL of 1X Assay Diluent into remaining tubes.

10.6 Prepare **Standard #2** by adding 120 µL Standard #1 to tube #2 and mix thoroughly.

10.7 Prepare **Standard #3** by adding 120 µL Standard #2 to tube #3 and mix thoroughly.

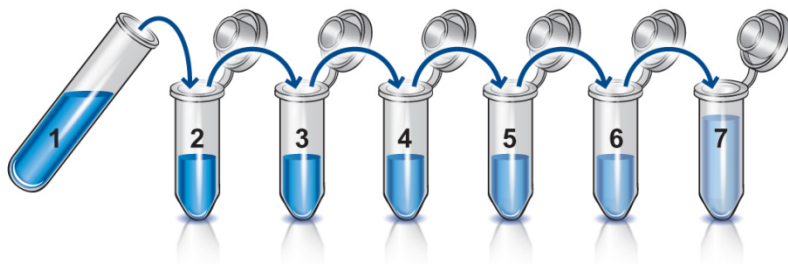
10.8 Using the table below as a guide, prepare further serial dilutions.

10.9 1X Assay Diluent 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	180	180	360	100,000	50,000
2	120	240	360	50,000	16,667
3	120	240	360	16,667	5,556
4	120	240	360	5,556	1,852
5	120	240	360	1,852	617.3
6	120	240	360	205.8	68.59
7	120	240	360	205.8	68.59
8	0	240	240	0	0



11. SAMPLE PREPARATION

General Sample Information:

- If your samples need to be diluted, Assay Diluent should be used for dilution of serum/plasma/culture supernatants
- Suggested dilution for normal serum/plasma: 10,000fold
- 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

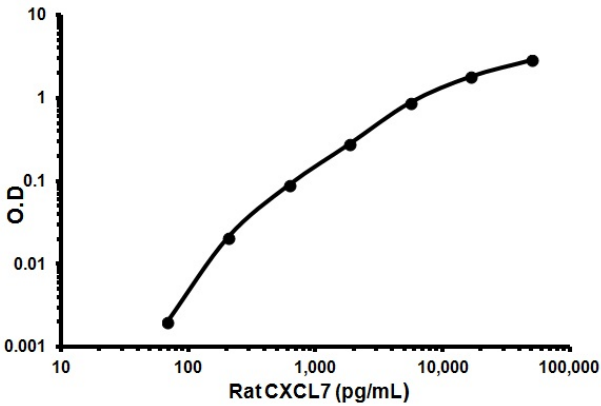
- **Equilibrate all materials and prepared reagents to room temperature (18 - 25°C) prior to use.**
 - **It is recommended to assay all standards, controls and samples in duplicate.**
- 13.1. Add 100 μ L of each standard and sample into appropriate wells. Cover well and incubate for 2.5 hours at room 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 μ L) using a multi-channel Pipette or auto washer. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
 - 13.3. Add 100 μ L of 1X Biotinylated CXCL7 Detection Antibody 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 μ L of 1X HRP-Streptavidin solution 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 μ L 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 μ L 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
68.59	0.002
205.8	0.021
617.3	0.089
1,852	0.279
5,556	0.881
16,667	1.818
50,000	2.869

16. TYPICAL SAMPLE VALUES

SENSITIVITY –

The minimum detectable dose of CXCL7 is typically less than 80 pg/mL.

RECOVERY –

Recovery was determined by spiking CXCL7 into normal Rat serum, plasma and cell culture media. Mean recoveries are as follows:

Sample Type	Average % Recovery	Range (%)
Serum	119.2	109-129
Plasma	116.3	93-132
Cell Culture Media	115.5	104-124

LINEARITY OF DILUTION -

Serum Dilution	Average % Expected Value	Range (%)
1:2	123.4	114-133
1:4	126.4	117-140

Plasma Dilution	Average % Expected Value	Range (%)
1:2	111.4	102-120
1:4	97.30	86-105

Cell Culture Media Dilution	Average % Expected Value	Range (%)
1:2	112.5	102-121
1:4	88.92	80-97

PRECISION –

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

17. ASSAY SPECIFICITY

Cross Reactivity: This ELISA kit shows no cross-reactivity with the following cytokines tested: rat CINC-2, CINC-3, CNTF, Fractalkine, IL-1a, IL-1b, IL-4, IL-6, IL-10, GM-CSF, IFN-g, Leptin, Lix, MCP-1, MIP-3a, b-NGF, TIMP-1, TNF-a.

18. 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.

19. 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