

Version 4a Last updated 15 March 2024

# ab219044 Rat CXCL1 ELISA Kit

For the quantitative measurement of rat CXCL1 in serum, plasma, cell culture supernatant, and cell extract samples.

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

# Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	5
9. Reagent Preparation	7
10. Standard Preparation	8
11. Sample Preparation	9
12. Plate Preparation	11
13. Assay Procedure	12
14. Calculations	14
15. Typical Data	15
16. Typical Sample Values	17
17. Assay Specificity	23
18. Species Reactivity	23
19. Troubleshooting	24
20. Notes	25

## 1. Overview

CXCL1 *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of rat CXCL1 protein in rat serum, plasma, cell culture supernatant, and cell extract samples.

The ELISA employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB Development Solution is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

CXCL1 (previously called GRO1 oncogene) is a member of the CXC family of chemokines. CXCL1 expression is reported in lymphocytes such as macrophages and neutrophils as well as epithelial cells. It plays a role in neutrophil chemotaxis. CXCL1 is involved in spinal cord development via inhibition of oligodendrocyte precursor migration as well angiogenesis, arteriogenesis, inflammation, wound healing and tumorigenesis.

Rat CXCL2 and CXCL3 share 65% and 64% sequence homology with rat CXCL1, respectively. Additionally, mouse CXCL1 is 89% homologous with rat CXCL1.

## 2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add 50  $\mu$ L standard or sample to appropriate wells



Add 50  $\mu$ L Capture Antibody Cocktail to all wells and incubate at room temperature for 30 minutes



Aspirate and wash each well three times with 350  $\mu$ L 1X Wash Buffer  
PT



Add 100  $\mu$ L Detector Antibody Cocktail to all wells and incubate at room temperature for 30 minutes



Aspirate and wash each well three times with 350  $\mu$ L 1X Wash Buffer  
PT



Add 100  $\mu$ L TMB Development Solution to each well and incubate for 10 minutes.



Add 100  $\mu$ L Stop Solution and read OD at 450 nm

### 3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

### 4. Storage and Stability

**Store kit at +4°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.**

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

## 5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

## 6. Materials Supplied

Item	Quantity	Storage Condition
Rat CXCL1 Capture Antibody 10X	600 µL	+4°C
Rat CXCL1 Detector Antibody 10X	600 µL	+4°C
Rat CXCL1 Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent CPI2	2 x 6 mL	+4°C
Wash Buffer PT 10X	2 x 20 mL	+4°C
Cell Extraction Buffer PTR 5X	10 mL	+4°C
Cell Extraction Enhancer Solution 50X	1 mL	+4°C
TMB Development Solution	12 mL	+4°C
Stop Solution	12 mL	+4°C
Sample Diluent NS	50 mL	+4°C
Anti-tag coated microplate (12 x 8 well strips)	96 Wells	+4°C
Plate Seal	1	+4°C

## 7. Materials Required, Not Supplied

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

- Microplate reader capable of measuring absorbance at 450 or 600 nm.
- Method for determining protein concentration (BCA assay recommended).
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for standard dilution.
- Plate shaker for all incubation steps.
- Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

## 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 is necessary to minimize background.
- As a guide, typical ranges of sample concentration for commonly used sample types are shown below in Sample Preparation (section 11).
- All samples should be mixed thoroughly and gently.
- Avoid multiple freeze/thaw of samples.
- Incubate ELISA plates on a plate shaker during all incubation steps.
- When generating positive control samples, it is advisable to change pipette tips after each step.
- The provided Cell Extraction Enhancer Solution 50X may precipitate when stored at + 4°C. To dissolve, warm briefly at + 37°C and mix gently. The Cell Extraction Enhancer Solution 50X can be stored at room temperature to avoid precipitation.
- **To avoid high background always add samples or standards to the well before the addition of the antibody cocktail.**
- **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. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.
- Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

### 9.1 1X Cell Extraction Buffer PTR (For cell and tissue extracts only):

Prepare 1X Cell Extraction Buffer PTR by diluting Cell Extraction Buffer PTR 5X and 50X Cell Extraction Enhancer Solution to 1X with deionized water. To make 10 mL 1X Cell Extraction Buffer PTR combine 7.8 mL deionized water, 2 mL Cell Extraction Buffer PTR 5X and 200  $\mu$ L Cell Extraction Enhancer Solution 50X. Mix thoroughly and gently. If required protease inhibitors can be added.

Alternative – Enhancer may be added to 1X Cell Extraction Buffer PTR after extraction of cells or tissue. Refer to note in the Troubleshooting section.

### 9.2 1X Wash Buffer PT:

Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 110 mL 1X Wash Buffer PT combine 11 mL Wash Buffer PT 10X with 99 mL deionized water. Mix thoroughly and gently.

### 9.3 Capture Antibody:

Prepare Capture Antibody by diluting the 10X Capture Antibody in Sample Diluent NS and Antibody Diluent CPI2. To make 3 mL of the Capture Antibody Cocktail combine 300  $\mu$ L 10X Capture Antibody and 300  $\mu$ L Sample Diluent NS with 2.4 mL Antibody Diluent CPI2. Mix thoroughly and gently.

### 9.4 Detector Antibody:

Prepare Detector Antibody by diluting the 10X Detector Antibody in Sample Diluent NS and Antibody Diluent CPI2. To make 6 mL of the Detector Antibody Cocktail combine 3.3 mL Sample Diluent NS and 300  $\mu$ L 10X Detector Antibody into 2.4 mL Antibody Diluent CPI2. Mix thoroughly and gently.

## 10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

**10.1 IMPORTANT:** If the protein standard vial has a volume identified on the label, reconstitute the rat CXCL1 protein standard by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the rat CXCL1 protein standard by adding 1,000  $\mu\text{L}$  Diluent. For **serum, plasma and cell culture supernatant sample measurements**, reconstitute the rat CXCL1 protein standard by adding Sample Diluent NS.

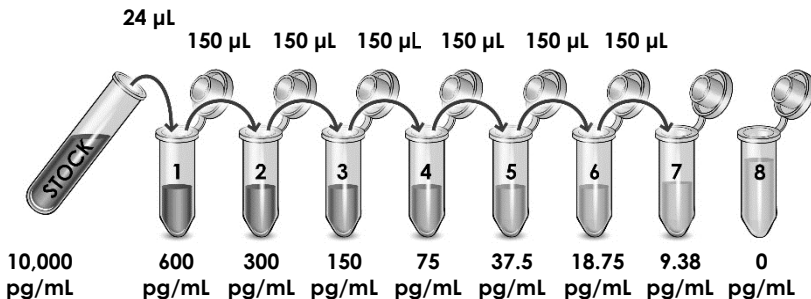
For **cell extract sample measurements**, reconstitute the rat CXCL1 protein standard by adding 1X Cell Extraction Buffer PTR.

Hold at room temperature for 10 minutes and mix thoroughly and gently. This is the 10,000  $\text{pg/mL}$  **Stock Standard** Solution.

**10.2** Label eight tubes, Standards 1–8.

**10.3** Add 376  $\mu\text{L}$  of appropriate diluent (see step 10.1) into tube number 1 and 150  $\mu\text{L}$  of appropriate diluent into numbers 2-8.

**10.4** Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



## 11. Sample Preparation

Typical Sample Dynamic Range	
Sample Type	Range
Plasma - Heparin	3.13 – 50%
Plasma - EDTA	3.13 – 50%
Plasma - Citrate	3.13 – 50%
Serum	3.13 – 50%
Lung Supernatant	1:1,600 – 1:100
Spleen Supernatant	1.56 – 25%
C6 Cell Extract	31.25 – 500 µg/mL

### 11.1 Plasma:

Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Dilute samples into Sample Diluent NS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

### 11.2 Serum:

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Dilute samples into Sample Diluent NS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

### 11.3 Cell Culture Supernatants:

Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants and dilute samples into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

## **11.4 Preparation of extracts from cell pellets:**

- 11.4.1 Collect non-adherent cells by centrifugation or scrape to collect adherent cells from the culture flask. Typical centrifugation conditions for cells are 500 x g for 5 minutes at 4°C.
- 11.4.2 Rinse cells twice with PBS.
- 11.4.3 Solubilize pellet at  $2 \times 10^7$  cell/mL in chilled 1X Cell Extraction Buffer PTR.
- 11.4.4 Incubate on ice for 20 minutes.
- 11.4.5 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.4.6 Transfer the supernatants into clean tubes and discard the pellets.
- 11.4.7 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.4.8 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

## **11.5 Preparation of extracts from adherent cells by direct lysis (alternative protocol):**

- 11.5.1 Remove growth media and rinse adherent cells 2 times in PBS.
- 11.5.2 Solubilize the cells by addition of chilled 1X Cell Extraction Buffer PTR directly to the plate (use 750  $\mu$ L - 1.5 mL 1X Cell Extraction Buffer PTR per confluent 15 cm diameter plate).
- 11.5.3 Scrape the cells into a microfuge tube and incubate the lysate on ice for 15 minutes.
- 11.5.4 Centrifuge at 18,000 x g for 20 minutes at 4°C.
- 11.5.5 Transfer the supernatants into clean tubes and discard the pellets.
- 11.5.6 Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay.
- 11.5.7 Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

## 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 plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.
- For each assay performed, a minimum of two wells must be used as the zero control.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Differences in well absorbance or “edge effects” have not been observed with this assay.

## 13. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
  - We recommend that you assay all standards, controls and samples in duplicate.
- 13.1** Prepare all reagents, working standards, and samples as directed in the previous sections.
  - 13.2** Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
  - 13.3** Add 50 µL of all sample or standard to appropriate wells.
  - 13.4** Add 50 µL of the Capture Antibody to each well.
  - 13.5** Seal the plate and incubate for 30 minutes at room temperature on a plate shaker set to 400 rpm.
  - 13.6** Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
  - 13.7** Add 100 µL of the Detector Antibody Cocktail to each well.
  - 13.8** Seal the plate and incubate for 30 minutes at room temperature on a plate shaker set to 400 rpm.
  - 13.9** Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
  - 13.10** Add 100 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm.

*Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes. Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the*

*high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.*

- 13.11** Add 100  $\mu$ L of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.
- 13.12** Alternative to 13.10 – 13.11: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

Mode	Kinetic
Wavelength:	600 nm
Time:	up to 20 min
Interval:	20 sec - 1 min
Shaking:	Shake between readings

**Δ Note:** that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100  $\mu$ L Stop Solution to each well and recording the OD at 450 nm.

- 13.13** Analyze the data as described below.

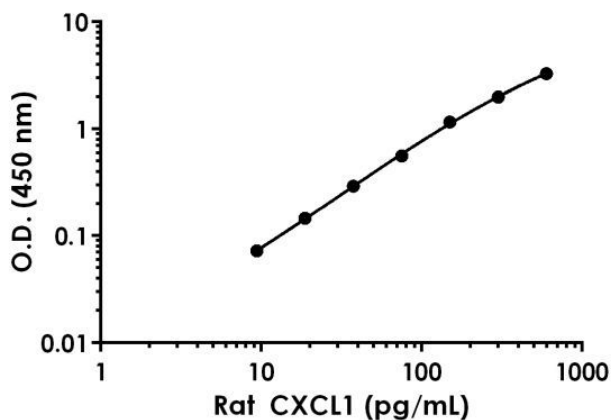
## 14. Calculations

- 14.1 Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
- 14.2 **Create a standard curve** by plotting the average blank control subtracted absorbance value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.
  - Δ **Note:** Most microplate reader software or graphing software will plot these values and fit a curve to the data. A four parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4 parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.
- 14.3 Determine the concentration of the target protein in the sample by interpolating the blank control subtracted **absorbance values against the standard curve**. Multiply the resulting value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.
- 14.4 Samples generating absorbance values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at an absorbance values less than that of the lowest standard should be retested in a less dilute form.



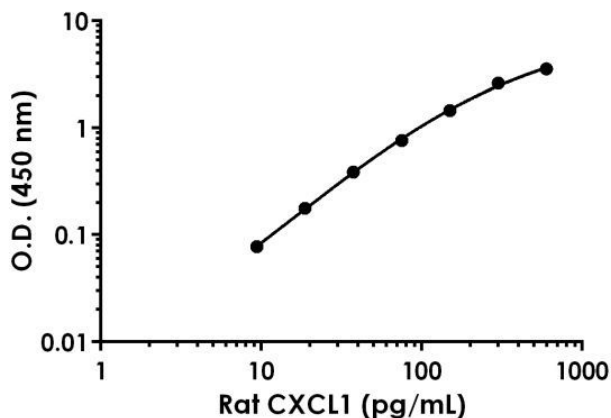
## 15. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.092	0.081	0.086
9.38	0.159	0.158	0.159
18.75	0.226	0.239	0.233
37.5	0.384	0.374	0.379
75	0.653	0.641	0.647
150	1.241	1.257	1.249
300	2.033	2.123	2.078
600	3.388	3.348	3.368

**Figure 1.** Example of rat CXCL1 standard curve in Sample Diluent NS. The CXCL1 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Standard Curve Measurements			
Concentration (pg/mL)	O.D 450 nm		Mean O.D
	1	2	
0	0.099	0.090	0.095
9.38	0.167	0.178	0.172
18.75	0.271	0.272	0.271
37.5	0.479	0.484	0.482
75	0.863	0.857	0.860
150	1.568	1.537	1.553
300	2.710	2.722	2.716
600	3.649	3.642	3.645

**Figure 2.** Example of rat CXCL1 standard curve in Sample Diluent 1X Cell Extraction Buffer PTR. The CXCL1 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

## 16. Typical Sample Values

### SENSITIVITY –

The calculated minimal detectable dose (MDD) was determined by calculating the mean of zero standard replicates and adding 2 standard deviations then extrapolating the corresponding concentration.

Sample Diluent Buffer	n=	Minimal Detectable Dose
Sample Diluent 10BS	24	2.13 pg/mL
1X Cell Extraction Buffer PTR	24	1.61 pg/mL

### RECOVERY –

Three concentrations of rat CXCL1 were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

Sample Type	Average % Recovery	Range (%)
Plasma – Heparin (50%)	114	110 – 118
Plasma – EDTA (50%)	106	103 – 108
Plasma – Citrate (50%)	97	88 – 112
Serum (50%)	99	88 – 109
Lung Supernatant (1:400)	101	85 – 116
Spleen Supernatant (10%)	96	90 – 100
RPMI Culture Media (25%)	95	90 – 102
C6 Cell Extract (125 µg/mL)	103	100 - 104

## Linearity of Dilution

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Native CXCL1 was measured in rat lung and spleen supernatant and rat C6 cell extract samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS for supernatant samples. Sample dilutions are made in Sample Diluent 1X Cell Extraction Buffer PTR for extract samples.

Recombinant rat CXCL1 was spiked into serum and plasma samples and diluted in a 2-fold dilution series in Sample Diluent NS.

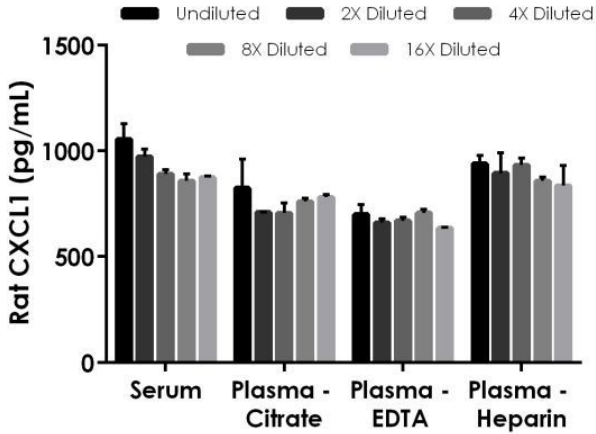
Dilution Factor	Interpolated value	50% Human Serum	50% Human Plasma (Citrate)	50% Human Plasma (EDTA)	50% Human Plasma (Heparin)
Undiluted	pg/mL	527.32	408.92	349.09	469.35
	<b>% Expected value</b>	100	100	100	100
2	pg/mL	243.19	177.19	165.19	222.68
	<b>% Expected value</b>	92	87	95	95
4	pg/mL	111.35	88.13	83.60	116.52
	<b>% Expected value</b>	84	86	96	99
8	pg/mL	53.63	47.55	44.17	53.64
	<b>% Expected value</b>	81	93	101	91
16	pg/mL	27.33	24.39	19.82	26.02
	<b>% Expected value</b>	83	95	91	89

Dilution Factor	Interpolated value	1% Lung Supernatant	25% Spleen Supernatant	500 µg/mL C6 Extract
Undiluted	pg/mL	559.46	239.48	525.06
	% Expected value	100	100	100
2	pg/mL	297.56	101.07	281.38
	% Expected value	106	84	107
4	pg/mL	153.35	48.79	127.94
	% Expected value	110	81	97
8	pg/mL	72.51	25.02	64.76
	% Expected value	104	84	99
16	pg/mL	33.54	11.92	34.00
	% Expected value	96	80	104

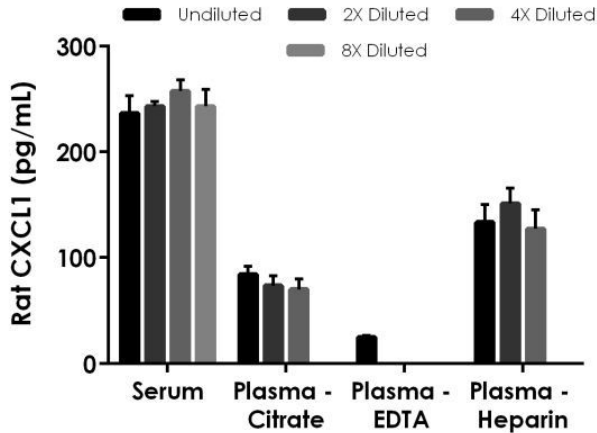
#### PRECISION –

Mean coefficient of variations of interpolated values from three concentrations of rat serum within the working range of the assay.

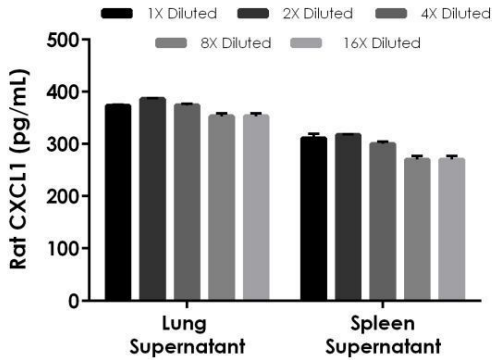
	Intra-Assay	Inter-Assay
n =	8	3
CV(%)	10.1	10.0



**Figure 3.** Interpolated concentrations of spiked CXCL1 in rat serum and plasma samples. The concentrations of CXCL1 were measured in duplicates, interpolated from the CXCL1 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (EDTA) 50%, and plasma (heparin) 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CXCL1 concentration was determined to be 930.66 pg/mL in neat serum, 756.22 pg/mL in neat plasma (citrate), 674.04 pg/mL in neat plasma (EDTA), and 891.86 pg/mL in neat plasma (heparin).

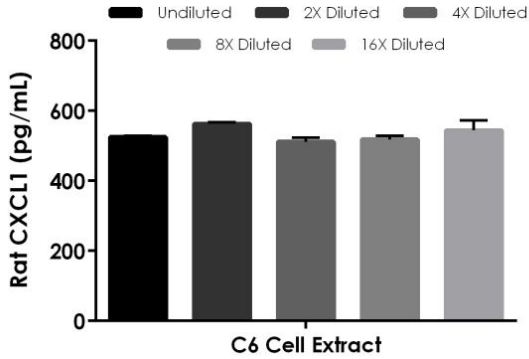


**Figure 4.** Interpolated concentrations of native CXCL1 in rat serum and plasma samples. The concentrations of CXCL1 were measured in duplicates, interpolated from the CXCL1 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (EDTA) 50%, and plasma (heparin) 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CXCL1 concentration was determined to be 245.16 pg/mL in neat serum, 75.79 pg/mL in neat plasma (citrate), 24.53 pg/mL in neat plasma (EDTA), and 137.19 pg/mL in neat plasma (heparin).



**Figure 5.** Interpolated concentrations of native CXCL1 in rat tissue culture supernatant samples. The concentrations of CXCL1 were measured in duplicates, interpolated from the CXCL1 standard curves and corrected for sample dilution. Undiluted samples are as follows: rat lung supernatant 1% and rat spleen supernatant 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CXCL1 concentration was determined to be 57,709.68 pg/mL in neat rat lung supernatant and 822.66 pg/mL in neat rat spleen supernatant. The rat lung supernatant was cultured for six days in RPMI base media with 10% Fetal Bovine Serum (FBS) plus 5 µg/mL of LPS. The rat spleen supernatant was cultured for three days in RPMI base media with 10% FBS.





**Figure 6.** Interpolated concentrations of native CXCL1 in rat C6 cell extract based on a 500  $\mu\text{g/mL}$  extract load. The concentrations of CXCL1 were measured in duplicate and interpolated from the CXCL1 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean  $\pm$  SD,  $n=2$ ). The mean CXCL1 concentration was determined to be 532.46  $\text{pg/mL}$  in C6 cell extract.

## 17. Assay Specificity

This kit recognizes both native and recombinant rat CXCL1 protein in serum, plasma, cell culture supernatant, and cell extract samples only.

### CROSS REACTIVITY

Recombinant rat CXCL2, mouse CXCL2, and mouse CXCL3 were prepared at 600 pg/mL and assayed for cross reactivity. No cross-reactivity was observed.

This kit is cross-reactive to mouse CXCL1 but is not suitable for use with mouse samples. Please contact our Technical Support team for more information.

### INTERFERENCE

Recombinant rat CXCL2 was prepared at 600 pg/mL and tested for interference. No interference with was observed.

## 18. Species Reactivity

This kit recognizes rat CXCL1 protein.

Other species reactivity was determined by measuring 1:2 diluted serum samples of various species, interpolating the protein concentrations from the rat standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in rat serum assayed at the same dilution.

Reactivity < 3% was determined for the following species:

- Human
- Cow

This kit is cross-reactive to mouse CXCL1 but is not suitable for use with mouse samples.

Please contact our Technical Support team for more information.

## 19. Troubleshooting

Problem	Reason	Solution
<b>Difficulty pipetting lysate; viscous lysate.</b>	Genomic DNA solubilized	Prepare 1X Cell Extraction Buffer PTR (without enhancer). Add enhancer to lysate after extraction.
<b>Poor standard curve</b>	Inaccurate Pipetting	Check pipettes
	Improper standard dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
<b>Low Signal</b>	Incubation times too brief	Ensure sufficient incubation times; increase to 2 or 3 hour standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
	Incubation times with TMB too brief	Ensure sufficient incubation time until blue color develops prior addition of Stop solution
<b>Large CV</b>	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
<b>Low sensitivity</b>	Improper storage of the ELISA kit	Store your reconstituted standards at -80°C, all other assay components 4°C. Keep TMB Development Solution protected from light.
<b>Precipitate in Diluent</b>	Precipitation and/or coagulation of components within the Diluent.	Precipitate can be removed by gently warming the Diluent to 37°C.

# Technical Support

Copyright © 2024 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)