



Cxcl16 (Mouse) ELISA Kit

Catalog Number KA0556

96 assays

Version: 07

Intended for research use only

www.abnova.com

Table of Contents

| | |
|---|-----------|
| Introduction | 3 |
| Intended Use | 3 |
| Background | 3 |
| Principle of the Assay | 3 |
| General Information | 4 |
| Materials Supplied | 4 |
| Storage Instruction | 4 |
| Materials Required but Not Supplied | 4 |
| Precautions for Use | 5 |
| Assay Protocol | 6 |
| Reagent Preparation | 6 |
| Sample Preparation | 7 |
| Assay Procedure | 7 |
| Data Analysis..... | 9 |
| Calculation of Results | 9 |
| Performance Characteristics | 9 |
| Resources | 10 |
| References | 10 |
| Plate Layout | 11 |

Introduction

Intended Use

For quantitative detection of mouse CXCL16 in cell culture supernates, serum and plasma (heparin, EDTA).

Background

Chemokine (C-X-C motif) ligand 16 (CXCL16) is a small cytokine belonging to the CXC chemokine family. Larger than other chemokines (with 254 amino acids), CXCL16 is composed of a CXC chemokine domain, a mucin-like stalk, a transmembrane domain and a cytoplasmic tail containing a potential tyrosine phosphorylation site that may bind SH2.¹ These are unusual features for a chemokine, and allow CXCL16 to be expressed as a cell surface bound molecule, as well as a soluble chemokine.² CXCL16 is produced by dendritic cells found in the T cell zones of lymphoid organs, and by cells found in the red pulp of the spleen.¹ Cells that bind and migrate in response to CXCL16 include several subsets of T cells, and natural killer T (NKT) cells.¹ CXCL16 interacts with the chemokine receptor CXCR6, also known as Bonzo.^{1,3} Expression of CXCL16 is induced by the inflammatory cytokines IFN-gamma and TNF-alpha.² The gene for human CXCL16 is located on chromosome 17p13.⁴ The standard product used in this kit is recombinant mouse CXCL16, consisting of 88 amino acids with the molecular mass of 9.9 KDa.

Principle of the Assay

The Cxcl16 (Mouse) ELISA kit was based on standard sandwich enzyme-linked immune-sorbent assay technology. A monoclonal antibody from rat specific for CXCL16 has been precoated onto 96-well plates. Standards and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for CXCL16 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the mouse CXCL16 amount of sample captured in plate.

General Information

Materials Supplied

List of component

| Component | Amount |
|---|-----------------|
| 96-well plate precoated with anti- mouse CXCL16 antibody. | 96 (8x12) wells |
| Lyophilized recombinant mouse CXCL16 standard | 10 ng/tube x 2 |
| Biotinylated anti- mouse CXCL16 antibody: dilution 1:100. | 130 μ L |
| Avidin-Biotin-Peroxidase Complex (ABC), dilution 1:100 | 130 μ L |
| Sample diluent buffer | 30 mL |
| Antibody diluent buffer | 12 mL |
| ABC diluent buffer | 12 mL |
| TMB color developing agent | 10 mL |
| TMB stop solution | 10 mL |

Storage Instruction

Store at 4°C for 6 months, at -20°C for 12 months. Avoid multiple freeze-thaw cycles.

Materials Required but Not Supplied

- ✓ Microplate reader in standard size.
- ✓ Automated plate washer.
- ✓ Adjustable pipettes and pipette tips. Multichannel pipettes are recommended in the condition of large amount of samples in the detection.
- ✓ Clean tubes and Eppendorf tubes.
- ✓ Washing buffer (neutral PBS or TBS).
- Preparation of 0.01 M TBS: Add 1.2 g Tris, 8.5 g NaCl; 450 μ L of purified acetic acid or 700 μ L of concentrated hydrochloric acid to 1000 mL H₂O and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1 L.
- Preparation of 0.01 M PBS: Add 8.5 g sodium chloride, 1.4 g Na₂HPO₄ and 0.2 g NaH₂PO₄ to 1000 mL distilled water and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1 L.

Precautions for Use

- ✓ To inspect the validity of experiment operation and the appropriateness of sample dilution proportion, pilot experiment using standards and a small number of samples is recommended.
- ✓ The TMB Color Developing agent is colorless and transparent before using, contact us freely if it is not the case.
- ✓ Before using the Kit, spin tubes and bring down all components to the bottom of tubes.
- ✓ Duplicate well assay is recommended for both standard and sample testing.
- ✓ Don't let 96-well plate dry, for dry plate will inactivate active components on plate.
- ✓ Don't reuse tips and tubes to avoid cross contamination.
- ✓ To avoid to use the reagents from different batches together.
- ✓ In order to avoid marginal effect of plate incubation due to temperature difference (reaction may be stronger in the marginal wells), it is suggested that the diluted ABC and TMB solution will be pre-warmed in 37°C for 30 min before using.

Assay Protocol

Reagent Preparation

- ✓ Reconstitution of the mouse CXCL16 standard: CXCL16 standard solution should be prepared no more than 2 hours prior to the experiment. Two tubes of CXCL16 standard (10 ng per tube) are included in each kit. Use one tube for each experiment.
 - 10,000 pg/mL of mouse CXCL16 standard solution: Add 1 mL sample diluent buffer into one tube, keep the tube at room temperature for 10 min and mix thoroughly.
 - 1000 pg/mL of mouse CXCL16 standard solution: Add 0.1 mL of the above 10 ng/mL CXCL16 standard solution into 0.9 mL sample diluent buffer and mix thoroughly.
 - 500 pg/mL → 15.6 pg/mL of mouse CXCL16 standard solutions: Label 6 Eppendorf tubes with 500 pg/mL, 250 pg/mL, 125 pg/mL, 62.5 pg/mL, 31.3 pg/mL, 15.6 pg/mL respectively. Aliquot 0.3 mL of the sample diluent buffer into each tube. Add 0.3 mL of the above 1000 pg/mL CXCL16 standard solution into 1st tube and mix. Transfer 0.3 mL from 1st tube to 2nd tube and mix. Transfer 0.3 mL from 2nd tube to 3rd tube and mix, and so on.

Note: The standard solutions are best used within 2 hours. The 10 ng/mL standard solution should be stored at 4°C for up to 12 hours, or at -20°C for up to 48 hours. Avoid repeated freeze-thaw cycles.

- ✓ Preparation of biotinylated anti-mouse CXCL16 antibody working solution: The solution should be prepared no more than 2 hours prior to the experiment.
 - The total volume should be: 0.1 mL/well x (the number of wells). (Allowing 0.1-0.2 mL more than total volume)
 - Biotinylated anti-mouse CXCL16 antibody should be diluted in 1:100 with the antibody diluent buffer and mixed thoroughly. (i.e. Add 1 µL Biotinylated anti-mouse CXCL16 antibody to 99 µL antibody diluent)

- ✓ Preparation of Avidin-Biotin-Peroxidase Complex (ABC) working solution: The solution should be prepared no more than 1 hour prior to the experiment.
 - The total volume should be: 0.1 mL/well x (the number of wells). (Allowing 0.1-0.2 mL more than total volume)
 - Avidin- Biotin-Peroxidase Complex (ABC) should be diluted in 1:100 with the ABC dilution buffer and mixed thoroughly. (i.e. Add 1 µL ABC to 99 µL ABC diluent buffer.)

Sample Preparation

✓ Sample Preparation and Storage

Store samples to be assayed within 24 hours at 2-8°C. For long-term storage, aliquot and freeze samples at -20°C. Avoid repeated freeze-thaw cycles.

- Cell culture supernates: Remove particulates by centrifugation, assay immediately or aliquot and store samples at -20°C.
- Serum: Allow the serum to clot in a serum separator tube (about 4 hours) at room temperature. Centrifuge at approximately 1000 x g for 15 min. Analyze the serum immediately or aliquot and store samples at -20°C.
- Plasma: Collect plasma using heparin or EDTA as an anticoagulant. Centrifuge for 15 min at 1500 x g within 30 min of collection. Assay immediately or aliquot and store samples at -20°C.

✓ Sample Dilution Guideline

The user needs to estimate the concentration of the target protein in the sample and select a proper dilution factor so that the diluted target protein concentration falls near the middle of the linear regime in the standard curve. Dilute the sample using the provided diluent buffer. The following is a guideline for sample dilution. Several trials may be necessary in practice. The sample must be well mixed with the diluents buffer.

- High target protein concentration (10-100 ng/mL). The working dilution is 1:100. i.e. Add 1 µL sample into 99 µL sample diluent buffer.
- Medium target protein concentration (1-10 ng/mL). The working dilution is 1:10. i.e. Add 10 µL sample into 90 µL sample diluent buffer.
- Low target protein concentration (15.6-1000 pg/mL). The working dilution is 1:2. i.e. Add 50 µL sample to 50 µL sample diluent buffer.
- Very Low target protein concentration (≤ 15.6 pg/mL). No dilution necessary, or the working dilution is 1:2.

Assay Procedure

The ABC working solution and TMB color developing agent must be kept warm at 37°C for 30 min before use. When diluting samples and reagents, they must be mixed completely and evenly. Standard CXCL16 detection curve should be prepared for each experiment. The user will decide sample dilution fold by crude estimation of CXCL16 amount in samples.

1. Aliquot 0.1 mL per well of the 1000 pg/mL, 500 pg/mL, 250 pg/mL, 125 pg/mL, 62.5 pg/mL, 31.3 pg/mL, 15.6 pg/mL mouse CXCL16 standard solutions into the precoated 96-well plate. Add 0.1 mL of the sample diluent buffer into the control well (Zero well). Add 0.1 mL of each properly diluted sample of mouse cell culture supernates, serum or plasma (heparin, EDTA) to each empty well. See "Sample Dilution Guideline" above for details. It is recommended that each mouse CXCL16 standard solution and each sample be measured in duplicate.

2. Seal the plate with the cover and incubate at 37°C for 90 min.
3. Remove the cover, discard plate content, and blot the plate onto paper towels or other absorbent material. Do NOT let the wells completely dry at any time.
4. Add 0.1 mL of biotinylated anti-mouse CXCL16 antibody working solution into each well and incubate the plate at 37°C for 60 min.
5. Wash the plate 3 times with 0.01 M TBS or 0.01 M PBS, and each time let washing buffer stay in the wells for 1 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material. (Plate Washing Method: Discard the solution in the plate without touching the side walls. Blot the plate onto paper towels or other absorbent material. Soak each well with at least 0.3 mL PBS or TBS buffer for 1~2 minutes. Repeat this process two additional times for a total of THREE washes. *Note: For automated washing, aspirate all wells and wash THREE times with PBS or TBS buffer, overfilling wells with PBS or TBS buffer. Blot the plate onto paper towels or other absorbent material.*)
6. Add 0.1 mL of prepared ABC working solution into each well and incubate the plate at 37°C for 30 min.
7. Wash plate 5 times with 0.01 M TBS or 0.01 M PBS, and each time let washing buffer stay in the wells for 1-2 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material. (See Step 5 for plate washing method).
8. Add 90 µL of prepared TMB color developing agent into each well and incubate plate at 37°C in dark for 15-20 min (*Note: For reference only, the optimal incubation time should be determined by end user. And the shades of blue can be seen in the wells with the four most concentrated mouse CXCL16 standard solutions; the other wells show no obvious color.*).
9. Add 0.1 mL of prepared TMB stop solution into each well. The color changes into yellow immediately.
10. Read the O.D. absorbance at 450 nm in a microplate reader within 30 min after adding the stop solution.

✓ Summary

1. Add samples and standards and incubate the plate at 37°C for 90 min. Do not wash.
2. Add biotinylated antibodies and incubate the plate at 37°C for 60 min. Wash plate 3 times with 0.01 M TBS.
3. Add ABC working solution and incubate the plate at 37°C for 30 min. Wash plate 5 times with 0.01 M TBS.
4. Add TMB color developing agent and incubate the plate at 37°C in dark for 15-20 min.
5. Add TMB stop solution and read.

Data Analysis

Calculation of Results

For calculation, (the relative O.D.₄₅₀) = (the O.D.₄₅₀ of each well) – (the O.D.₄₅₀ of Zero well). The standard curve can be plotted as the relative O.D.₄₅₀ of each standard solution (Y) vs. the respective concentration of the standard solution (X). The mouse CXCL16 concentration of the samples can be interpolated from the standard curve.

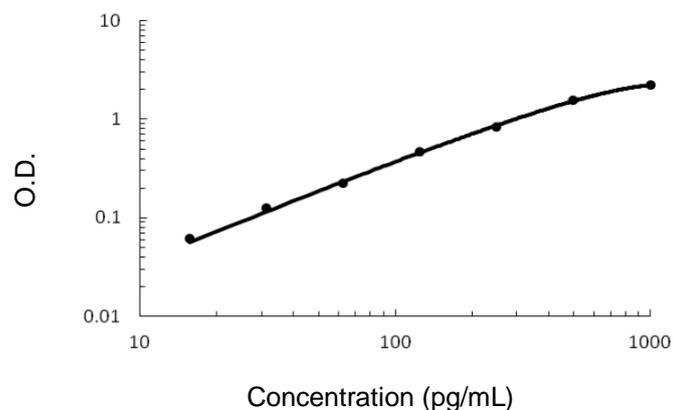
Note: if the samples measured were diluted, multiply the dilution factor to the concentrations from interpolation to obtain the concentration before dilution.

✓ Typical result

Typical Data Obtained from mouse CXCL16

(TMB reaction incubate at 37°C for 17 min)

| | | | | | | | | |
|-----------------------|-------|-------|-------|-------|-------|-------|-------|-------|
| Concentration (pg/mL) | 0.0 | 15.6 | 31.2 | 62.5 | 125 | 250 | 500 | 1000 |
| O.D | 0.013 | 0.062 | 0.125 | 0.223 | 0.465 | 0.841 | 1.547 | 2.216 |



This standard curve was generated for demonstration purpose only. A standard curve must be run with each assay.

Performance Characteristics

- ✓ Range: 15.6 pg/mL – 1000 pg/mL
- ✓ Sensitivity: < 1 pg/mL
- ✓ Specificity: Natural and recombinant mouse CXCL16
- ✓ Cross-reactivity: No detectable cross-reactivity with other relevant proteins

Resources

References

1. Matloubian M, David A, Engel S, Ryan J, Cyster J (2000). "A transmembrane CXC chemokine is a ligand for HIV-coreceptor Bonzo". *Nat Immunol* 1 (4): 298–304.
2. Abel S, Hundhausen C, Mentlein R, Schulte A, Berkhout T, Broadway N, Hartmann D, Sedlacek R, Dietrich S, Muetze B, Schuster B, Kallen K, Saftig P, Rose-John S, Ludwig A (2004). "The transmembrane CXC-chemokine ligand 16 is induced by IFN-gamma and TNF-alpha and shed by the activity of the disintegrin-like metalloproteinase ADAM10". *J Immunol* 172 (10): 6362–72.
3. Wilbanks A, Zondlo S, Murphy K, Mak S, Soler D, Langdon P, Andrew D, Wu L, Briskin M (2001). "Expression cloning of the STRL33/BONZO/TYMSTR ligand reveals elements of CC, CXC, and CX3C chemokines". *J Immunol* 166 (8): 5145–54.
4. Matloubian, M.; David, A.; Engel, S.; Ryan, J. E.; Cyster, J. G. : A transmembrane CXC chemokine is a ligand for HIV-coreceptor Bonzo. *Nature Immun.* 1: 298-304, 2000.

Plate Layout

| | | | | | | | | |
|----|---|---|---|---|---|---|---|---|
| 12 | | | | | | | | |
| 11 | | | | | | | | |
| 10 | | | | | | | | |
| 9 | | | | | | | | |
| 8 | | | | | | | | |
| 7 | | | | | | | | |
| 6 | | | | | | | | |
| 5 | | | | | | | | |
| 4 | | | | | | | | |
| 3 | | | | | | | | |
| 2 | | | | | | | | |
| 1 | | | | | | | | |
| | A | B | C | D | E | F | G | H |