



BSG (Human) ELISA Kit

Catalog Number KA0988

96 assays

Version: 04

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 | 6 |
| 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 human Emmprin in cell culture supernates, serum, plasma(heparin, EDTA), saliva, urine and human milk.

Background

Basigin is a member of the immunoglobulin superfamily that is also known as Emmprin, short for extracellular matrix metalloproteinase inducer, and recently has been designated CD147 (cluster of differentiation 147).^{[1][2]} It is a member of the immunoglobulin superfamily, with a structure related to the putative primordial form of the family. As members of the immunoglobulin superfamily play fundamental roles in intercellular recognition involved in various immunologic phenomena, differentiation, and development, basigin is thought also to play a role in intercellular recognition (Miyauchi et al., 1991; Kanekura et al., 1991). This protein is a determinant for the Ok blood group system. Basigin is a type I integral membrane receptor that has many ligands, including the cyclophilin (CyP) proteins Cyp-A and CyP-B and certain integrins.^{[3][4][5]} It is expressed by many cell types, including epithelial cells, endothelial cells and leukocytes. The standard product used in this kit is recombinant human Emmprin · A22-H205 - 47.4KDa.

Principle of the Assay

BSG (Human) ELISA Kit was based on standard sandwich enzyme-linked immune-sorbent assay technology. Human Emmprin specific-specific polyclonal antibodies were precoated onto 96-well plates. The human specific detection monoclonal antibodies were biotinylated. The test samples and biotinylated detection antibodies were added to the wells subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to 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 human BSG amount of sample captured in plate.

General Information

Materials Supplied

List of component

| Component | Amount |
|---|-----------|
| 96-well plate precoated with anti- human Emmprin antibody | 1 plate |
| Lyophilized recombinant human Emmprin standard | 10 ng × 2 |
| Biotinylated anti- human Emmprin antibody, dilution 1:100 | 130 µl |
| Avidin-Biotin-Peroxidase Complex (ABC), dilution 1:100 | 130 µl |
| 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 frequent use, at -20°C for infrequent use. 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 0.01M TBS: Add 1.2g Tris, 8.5g NaCl; 450µl of purified acetic acid or 700µl of concentrated hydrochloric acid to 1000ml H₂O and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

Preparation of 0.01 M PBS: Add 8.5g sodium chloride, 1.4g Na₂HPO₄ and 0.2g NaH₂PO₄ to 1000ml distilled water and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

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 human Emmprin standard: Emmprin standard solution should be prepared no more than 2 hours prior to the experiment. Two tubes of Emmprin standard (10ng per tube) are included in each kit. Use one tube for each experiment.
 - 10,000pg/ml of human Emmprin standard solution: Add 1 ml sample diluent buffer into one tube, keep the tube at room temperature for 10 min and mix thoroughly.
 - 2000pg/ml of human Emmprin standard solution: Add 0.2 ml of the above 10ng/ml Emmprin standard solution into 0.8 ml sample diluent buffer and mix thoroughly.
 - 1000pg/ml→31.2pg/ml of human Emmprin standard solutions: Label 6 Eppendorf tubes with 1000pg/ml, 500pg/ml, 250pg/ml, 125pg/ml, 62.5pg/ml, 31.2pg/ml, respectively. Aliquot 0.3 ml of the sample diluent buffer into each tube. Add 0.3 ml of the above 2000pg/ml Emmprin 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 may 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-human Emmprin antibody working solution: The solution should be prepared no more than 2 hours prior to the experiment.
 - The total volume should be: 0.1ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume)
 - Biotinylated anti-human Emmprin antibody should be diluted in 1:100 with the antibody diluent buffer and mixed thoroughly. (i.e. Add 1µl Biotinylated anti-human Emmprin antibody to 99µl antibody diluent buffer.)
- ✓ 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: 0.1ml/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 1µl ABC to 99µl 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.
- Saliva: Collect saliva using a collection device without any protein binding or filtering capabilities such as a Salivette or aliquot and store samples at -20°C.
- Urine: Aseptically collect the first urine of the day, micturate directly into a sterile container. Remove particular impurities by centrifugation, assay immediately or aliquot and store samples at -20°C.
- Human milk: Centrifuge for 15 min at 1500 x g at 2-8°C. Collect the aqueous fraction and repeat this process 3 times. Filter through a 0.2µm filter and assay immediately or aliquot and store samples at -80°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 (20-200ng/ml). The working dilution is 1:100. i.e. Add 1 µl sample into 99 µl sample diluent buffer.
- Medium target protein concentration (2-20ng/ml). The working dilution is 1:10. i.e. Add 10 µl sample into 90 µl sample diluent buffer.
- Low target protein concentration (31.2-2000pg/ml). The working dilution is 1:2. i.e. Add 50 µl sample to 100 µl sample diluent buffer.
- Very Low target protein concentration (≤ 31.2 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 Emmprin detection curve should be prepared for each experiment. The user will decide sample dilution fold by crude estimation of Emmprin amount in samples.

1. Aliquot 0.1ml per well of the 2000pg/ml, 1000pg/ml, 500pg/ml, 250pg/ml, 125pg/ml, 62.5pg/ml, 31.2pg/ml human Emmprin standard solutions into the precoated 96-well plate. Add 0.1ml of the sample diluent buffer into the control well (Zero well). Add 0.1ml of each properly diluted sample of human cell culture

supernates, serum, plasma(heparin, EDTA), saliva, urine or human milk to each empty well. See “Sample Dilution Guideline” above for details. It is recommend that each human Emmprin standard solution and each sample is 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 0.1ml of biotinylated anti-human Emmprin antibody working solution into each well and incubate the plate at 37°C for 60 min.
5. Wash the plate 3 times with 0.01M TBS or 0.01M 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 0.1ml of prepared ABC working solution into each well and incubate the plate at 37°C for 30 min.
7. Wash plate 5 times with 0.01M TBS or 0.01M 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 human Emmprin standard solutions; the other wells show no obvious color*).
9. Add 0.1ml of prepared TMB stop solution into each well. The color changes into yellow immediately.
10. Read the O.D. absorbance at 450nm 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 0.01M TBS.
3. Add ABC working solution and incubate the plate at 37°C for 30 min. Wash plate 5 times with 0.01M 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 human Emmprin 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 Human Emmprin

| Concentration pg/ml | 0.0 | 31.2 | 62.5 | 125 | 250 | 500 | 1000 | 2000 |
|------------------------|-------|-------|-------|-------|-------|-------|-------|-------|
| O.D. | 0.040 | 0.099 | 0.160 | 0.281 | 0.472 | 0.819 | 1.583 | 2.156 |

(TMB reaction incubate at 37°C for 15 min)

Performance Characteristics

- ✓ Range
31.2pg/ml-2000pg/ml
- ✓ Sensitivity
< 2pg/ml
- ✓ Specificity
Natural and recombinant human Emmprin.
- ✓ Cross-reactivity
No detectable cross-reactivity with other relevant proteins

Resources

References

1. Kasinrerker W, Fiebigler E, Stefanová I, Baumruker T, Knapp W, Stockinger H (1992). "Human leukocyte activation antigen M6, a member of the Ig superfamily, is the species homologue of rat OX-47, mouse basigin, and chicken HT7 molecule". *J Immunol* 149 (3): 847–54.
2. Yurchenko V, Constant S, Bukrinsky M (2006). "Dealing with the family: CD147 interactions with cyclophilins". *Immunology* 117 (3): 301–9.
3. Yurchenko V, Zybarth G, O'Connor M, Dai W, Franchin G, Hao T, Guo H, Hung H, Toole B, Gallay P, Sherry B, Bukrinsky M (2002). "Active site residues of cyclophilin A are crucial for its signaling activity via CD147". *J Biol Chem* 277 (25): 22959–65.
4. Yurchenko V, O'Connor M, Dai W, Guo H, Toole B, Sherry B, Bukrinsky M (2001). "CD147 is a signaling receptor for cyclophilin B". *Biochem Biophys Res Commun* 288 (4): 786–8.
5. Berditchevski F, Chang S, Bodorova J, Hemler M (1997). "Generation of monoclonal antibodies to integrin-associated proteins. Evidence that alpha3beta1 complexes with BSG/basigin/OX47/M6". *J Biol Chem* 272 (46): 29174–80.

Plate Layout

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