
Product Manual

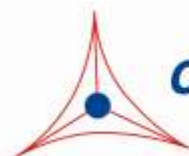
Homocysteine ELISA Kit

Catalog Number

STA-670

96 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Homocysteine is an amino acid intermediate formed during the production of the essential dietary amino acid methionine (Figure 1). Homocysteine is a homologue of cysteine, differing from cysteine only in that it contains an extra side chain methylene bridge. About 80% of homocysteine found in plasma is bound to protein. High levels of homocysteine in the blood have been associated with premature incidences of vascular disease, and homocysteine is likely to be a risk factor for heart disease. Homocysteine initially stimulates the production of nitric oxide in endothelial cells but ultimately reduces nitric oxide bioavailability and increases oxidative stress by blocking glutathione peroxidase activity as well as causing cellular oxidative degradation (increasing free radical generation). In addition, elevated homocysteine levels leads to increased platelet and leukocyte adhesion and activation, increased vasoconstriction, and increased proliferation of smooth muscle (a hallmark of atherosclerosis).

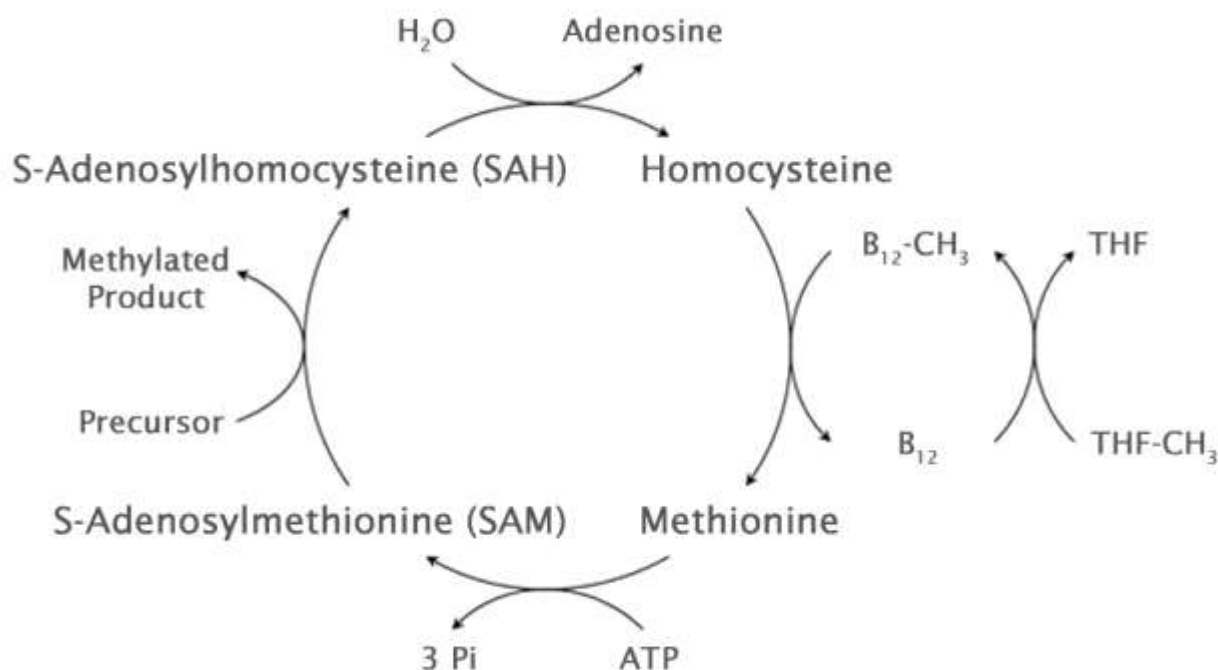


Figure 1. Metabolism of homocysteine.

Cell Biolabs' Homocysteine ELISA Kit is a competitive enzyme immunoassay developed for the detection and quantitation of homocysteine in plasma, serum, lysates, or other biological fluid samples. The kit has a detection sensitivity limit of 10 ng/mL Homocysteine-BSA. Each kit provides sufficient reagents to perform up to 96 assays including standard curve and unknown samples.

Related Products

1. STA-361: Human ApoAI and ApoB Duplex ELISA Kit
2. STA-369: OxiSelect™ Human Oxidized LDL ELISA Kit (MDA-LDL Quantitation)

3. STA-390: Total Cholesterol Assay Kit
4. STA-391: HDL and LDL/VLDL Cholesterol Assay Kit
5. STA-396: Serum Triglyceride Quantification Kit (Colorimetric)
6. STA-398: Free Glycerol Assay Kit (Colorimetric)
7. STA-671: S-Adenosylhomocysteine (SAM) ELISA Kit
8. STA-672: S-Adenosylmethionine (SAM) ELISA Kit

Kit Components

Box 1 (shipped at room temperature)

1. 96 Well Protein Binding Plate (Part No. 231001): One strip well 96 well plate.
2. Anti-Homocysteine Antibody (500X) (Part No. 267002): One 15 μ L vial.
3. Secondary Antibody, HRP Conjugate (1000X) (Part No. 231009): One 20 μ L vial.
4. Assay Diluent (Part No. 310804): One 50 mL bottle.
5. 10X Wash Buffer (Part No. 310806): One 100 mL bottle.
6. Substrate Solution (Part No. 310807): One 12 mL amber bottle.
7. Stop Solution (Part. No. 310808): One 12 mL bottle.

Box 2 (shipped on blue ice packs)

1. Homocysteine Conjugate (1000X) (Part No. 267001): One 20 μ L vial.
2. Homocysteine-BSA Standard (Part No. 267003): One 20 μ L vial of 4 mg/mL homocysteine conjugated to BSA in PBS.

Materials Not Supplied

1. Plasma, serum, or other biological fluids
2. Phosphate Buffered Saline (PBS)
3. PBS containing 0.1% Bovine Serum Albumin (BSA)
4. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
5. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
6. Multichannel micropipette reservoir
7. Microplate reader capable of reading at 450 nm (620 nm as optional reference wave length)

Storage

Upon receipt, store Homocysteine Conjugate (1000X) and Homocysteine-BSA Standard at -20°C. Store the rest of the kit at 4°C.

Preparation of Reagents

- Homocysteine Conjugate Coated Plate: Determine the number of wells to be used, and dilute the Homocysteine Conjugate 1:1000 into PBS. Add 100 µL of 1X homocysteine conjugate to each well of the 96-well Protein Binding Plate. Incubate for 2 hrs at 37°C or overnight at 4°C. Remove the diluted homocysteine conjugate, blotting plate on paper towels to remove excess fluid. Wash wells 3 times with 200 µL of PBS and blot on paper towels to remove excess fluid. Add 200 µL of Assay Diluent to each well and block for 1 hour at room temperature. Transfer the plate to 4°C until ready to begin the assay.

Note: The Homocysteine Conjugate Coated Plate is not stable long-term. We recommend using it within 24 hours after coating.

- 1X Wash Buffer: Dilute the 10X Wash Buffer to 1X with deionized water. Stir to homogeneity.
- Anti-Homocysteine Antibody and Secondary Antibody, HRP Conjugate: Immediately before use dilute the Anti-Homocysteine Antibody 1:500 and the Secondary Antibody, HRP Conjugate 1:1000 with Assay Diluent. Do not store diluted solutions.

Preparation of Standard Curve

Prepare a dilution series of Homocysteine-BSA standards in the concentration range of 0 to 40 µg/mL in Assay Diluent (Table 1).

Standard Tubes	4 mg/mL Homocysteine-BSA Standard (µL)	Assay Diluent (µL)	Homocysteine-BSA (µg/mL)
1	4	396	40
2	100 of Tube #1	300	10
3	100 of Tube #2	300	2.5
4	100 of Tube #3	300	0.625
5	100 of Tube #4	300	0.156
6	100 of Tube #5	300	0.039
7	100 of Tube #6	300	0.010
8	0	300	0

Table 1. Preparation of Homocysteine-BSA Standards.

Preparation of Samples

The following recommendations are only guidelines and may be altered to optimize or complement the user's experimental design.

- Plasma: Collect blood with heparin or EDTA and centrifuge for 10 minutes at 1000 g at 4°C. Remove the plasma and assay immediately or store samples at -80°C for up to three months. Normal plasma samples should be diluted 2- to 10-fold with PBS containing 0.1% BSA immediately before running the ELISA.

- Serum: Harvest serum and centrifuge for 10 minutes at 1000 g at 4°C. Assay immediately or store samples at -80°C for up to three months. Normal serum samples should be diluted 2- to 10-fold with PBS containing 0.1% BSA immediately before running the ELISA.
- Tissue homogenate: Weigh and homogenize the tissue on ice in 5-10 mL cold PBS per gram of tissue. Centrifuge at 10,000 x g for 15 minutes at 4°C. Remove the supernatant and store on ice. Store any unused supernatant at -80°C for up to three months.
- Cell lysate: Collect cells by centrifuging at 2000 x g for 10 minutes at 4°C. Sonicate or homogenize the cell pellet on ice in 1-2 mL cold PBS. Centrifuge at 10,000 x g for 15 minutes at 4°C. Remove the supernatant and store on ice. Aliquot and store the supernatant for use in the assay. Store any unused supernatant at -80°C for up to three months.
- Other biological fluids: Centrifuge samples for 10 minutes at 1000 g at 4°C and recover supernatant. Assay immediately or store samples at -80°C for up to three months.

Assay Protocol

1. Prepare and mix all reagents thoroughly before use.
2. Each unknown sample (see Preparation of Samples section), Homocysteine-BSA standard, and blank should be assayed in duplicate.
3. Remove the Assay Diluent from the plate and add 50 µL of unknown sample or standard to the Homocysteine Conjugate Coated Plate. Incubate at room temperature for 10 minutes on an orbital shaker.
4. Add 50 µL of diluted Anti-Homocysteine Antibody (see Preparation of Reagents section) to each well. Incubate at room temperature for 1 hour on an orbital shaker.
5. Wash microwell strips 3 times with 250 µL 1X Wash Buffer per well with thorough aspiration between each wash. After each wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
6. Add 100 µL of the diluted Secondary Antibody, HRP Conjugate to each well. Incubate at room temperature for 1 hour on an orbital shaker. During this incubation, warm Substrate Solution to room temperature.
7. Wash the strip wells 3 times according to step 5 above. Proceed immediately to the next step.
8. Add 100 µL of Substrate Solution to each well, including the blank wells. Incubate at room temperature on an orbital shaker. Actual incubation time may vary from 2-30 minutes.

Note: Watch plate carefully; if color changes rapidly, the reaction may need to be stopped sooner to prevent saturation.

9. Stop the enzyme reaction by adding 100 μL of Stop Solution into each well, including the blank wells. Results should be read immediately (color will fade over time).
10. Read absorbance of each microwell on a spectrophotometer using 450 nm as the primary wave length.

Example of Results

The following figures demonstrate typical Homocysteine Competitive ELISA Kit results. One should use the data below for reference only. This data should not be used to interpret actual results.

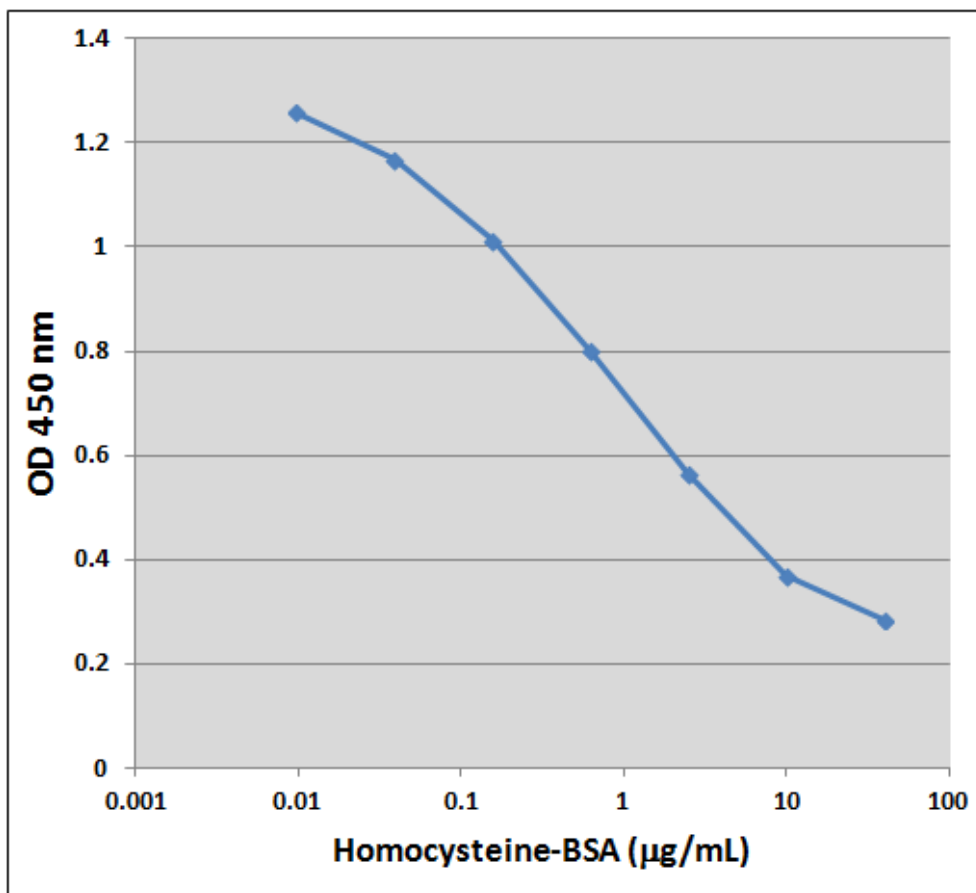


Figure 2: Homocysteine-BSA Standard Curve.

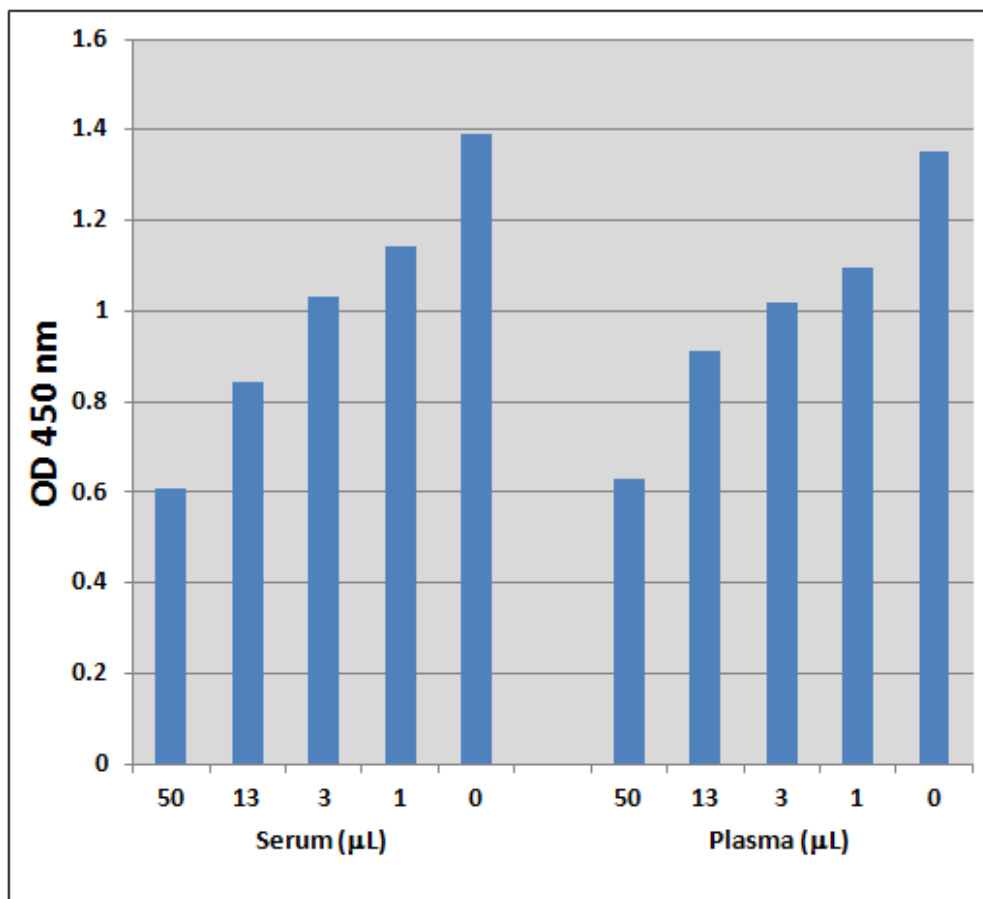


Figure 3: Homocysteine Detection in Human Serum and Plasma.

References

1. Thambyrajah J, and Townend JN (2000) *European Heart J.* **21**:967-974.
2. Starkebaum G and Harlan JM. (1986) *J. Clin Invest.* **77**:1370-6.
3. Loscalzo J (1996) *J. Clin Invest.* **98**:5-7.
4. Upchurch GR, Jr., Welch GN Fabian AJ, Freedman JE, Johnson JL, Keaney JF Jr, and Loscalzo J. (1997) *J. Biol Chem* **272**:17012-17017.
5. Tyagi SC (1998) *Am. J. Physiol.* **274** (2 Pt 1): C396-405.

Recent Product Citations

1. Ali, M.M. et al. (2021). DNA methylation profile of genes involved in inflammation and autoimmunity correlates with vascular function in morbidly obese adults. *Epigenetics*. doi: 10.1080/15592294.2021.1876285.
2. Toriumi, K. et al. (2021). Combined glyoxalase 1 dysfunction and vitamin B6 deficiency in a schizophrenia model system causes mitochondrial dysfunction in the prefrontal cortex. *Redox Biol.* **45**:102057. doi: 10.1016/j.redox.2021.102057.

3. Spechenkova, N. et al. (2021). The Resistance Responses of Potato Plants to Potato Virus Y Are Associated with an Increased Cellular Methionine Content and an Altered SAM:SAH Methylation Index. *Viruses*. **13**(6):955. doi: 10.3390/v13060955.
4. Cheng, C.K. et al. (2021). A high methionine and low folate diet alters glucose homeostasis and gut microbiome. *Biochem Biophys Res Commun*. doi: 10.1016/j.bbrep.2021.100921.
5. Fesenko, I. et al. (2020). Role of the methionine cycle in the temperature-sensitive responses of potato plants to potato virus Y. *Mol Plant Pathol*. doi: 10.1111/mpp.13009.
6. Siervo, M. et al. (2020). Nitrate-Rich Beetroot Juice Reduces Blood Pressure in Tanzanian Adults with Elevated Blood Pressure: A Double-Blind Randomized Controlled Feasibility Trial. *J Nutr*. doi: 10.1093/jn/nxaa170.
7. Alfarisi, H.A.H. et al. (2020). Anti-inflammatory Effects of Trihoney in Hypercholesterolemic Atherosclerotic Rabbits: A Comparative Study With Atorvastatin. *Mal J Med Health Sci*. **16**(2): 230-236.
8. Fairbairn, P. et al. (2020). Effects of a high DHA multi-nutrient supplement and exercise on mobility and cognition in older women (MOBILE): A randomised semi-blinded placebo controlled study. *Br J Nutr*. doi: 10.1017/S0007114520000719.
9. Chen, Y.C. et al. (2019). Melatonin rescued methotrexate-induced spatial deficit and hyperhomocysteinemia and increased asymmetric dimethylarginine in plasma and dorsal hippocampus in developing rats. *Life Sci*. doi: 10.1016/j.lfs.2019.116931.
10. Fan, Y. et al. (2019). Excessive Methionine Supplementation Exacerbates the Development of Abdominal Aortic Aneurysm in Rats. *J Vasc Res*. doi: 10.1159/000501313.
11. Kang, Y. et al. (2019). Unraveling the role of quorum sensing-dependent metabolic homeostasis of the activated methyl cycle in a cooperative population of *Burkholderia glumae*. *Sci Rep*. **9**(1):11038. doi: 10.1038/s41598-019-47460-6.
12. Fairbairn, P. et al. (2019). Combining a high DHA multi-nutrient supplement with aerobic exercise: Protocol for a randomised controlled study assessing mobility and cognitive function in older women. *Prostaglandins Leukot Essent Fatty Acids*. **143**:21-30. doi: 10.1016/j.plefa.2019.04.001.
13. Campbell, M.S. et al. (2019). Influence of enhanced bioavailable curcumin on obesity-associated cardiovascular disease risk factors and arterial function: A double-blinded, randomized, controlled trial. *Nutrition*. **62**:135-139. doi: 10.1016/j.nut.2019.01.002.
14. May, J.L. et al. (2019). IDH3 α regulates one-carbon metabolism in glioblastoma. *Sci Adv*. **5**(1):eaat0456. doi: 10.1126/sciadv.aat0456.
15. Saadeh, N. et al. (2018). Serum homocysteine is associated with polycystic ovarian syndrome in Jordan. *Biomed Rep*. **9**(5):439-445. doi: 10.3892/br.2018.1149.
16. Zhou X. et al. (2018). Serine prevented high-fat diet-induced oxidative stress by activating AMPK and epigenetically modulating the expression of glutathione synthesis-related genes. *Biochim Biophys Acta Mol Basis Dis*. **1864**(2):488-498. doi: 10.1016/j.bbadis.2017.11.009.
17. Shah, T. et al. (2018). Vitamin B12 insufficiency in excess folic acid downregulates methylenetetrahydrofolate reductase gene and increases homocysteine, tumor necrosis factor-alpha, and oxidative stress in hepatocytes. *Indian J Health Sci Biomed Res*. **11**:12-8. doi: 10.4103/kleuhsj.kleuhsj_128_17.
18. Cao, G. et al. (2018). Effects of maternal serine supplementation on high-fat diet-induced oxidative stress and epigenetic changes in promoters of glutathione synthesis-related genes in offspring. *Journal of Functional Foods*. **47**:316-324. doi: 10.1016/j.jff.2018.05.067.

19. Toghil, B.J. et al. (2018). SMYD2 promoter DNA methylation is associated with abdominal aortic aneurysm (AAA) and SMYD2 expression in vascular smooth muscle cells. *Clin Epigenetics*. **10**:29. doi: 10.1186/s13148-018-0460-9.
20. Zhou, X. et al. (2018). Serine prevented high-fat diet-induced oxidative stress by activating AMPK and epigenetically modulating the expression of glutathione synthesis-related genes. *Biochim Biophys Acta*. **1864**(2):488-498. doi: 10.1016/j.bbadis.2017.11.009
21. Elamir, A. et al (2017). Role of long non-coding RNA CDKN2B-AS1 and Homocysteine in patients with ischemic cardiomyopathy. *WJPLS*. **3** (7): 11-18.
22. Shah, T. et al. (2016). Molecular and cellular effects of vitamin B12 forms on human trophoblast cells in presence of excessive folate. *Biomed Pharmacother*. **84**:526-534.
23. Liu, B. et al. (2016). A novel rat model of heart failure induced by high methionine diet showing evidence of association between hyperhomocysteinemia and activation of NF-kappaB. *Am J Transl Res*. **8**:117-124.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2015-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.