
Product Manual

OxiSelect™ HNE Adduct Competitive ELISA Kit, Trial Size

Catalog Number

STA-838-T

32 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Lipid peroxidation is a well-defined mechanism of cellular damage in animals and plants. Lipid peroxides are unstable indicators of oxidative stress in cells that decompose to form more complex and reactive compounds such as Malondialdehyde (MDA) and 4-hydroxynonenal (4-HNE), natural bi-products of lipid peroxidation. Oxidative modification of lipids can be induced *in vitro* by a wide array of pro-oxidant agents and occurs *in vivo* during aging and in certain disease conditions. Measuring the end products of lipid peroxidation is one of the most widely accepted assays for oxidative damage. These aldehydic secondary products of lipid peroxidation are generally accepted markers of oxidative stress.

Both MDA and HNE have been shown to be capable of binding to proteins and forming stable adducts, also termed advanced lipid peroxidation end products. These modifications of proteins by MDA or HNE can cause both structural and functional changes of oxidized proteins. Specifically, 4-HNE can react with lysine, histidine or cysteine residues in protein to form adducts.

Cell Biolabs' OxiSelect™ HNE Adduct Competitive ELISA Kit is an enzyme immunoassay developed for rapid detection and quantitation of HNE protein adducts. The quantity of HNE adduct in protein samples is determined by comparing its absorbance with that of a known HNE-BSA standard curve. Each Trial Size HNE Adduct Competitive ELISA Kit provides sufficient reagents to perform up to 32 assays, including standard curve and unknown protein samples.

Assay Principle

First, an HNE conjugate is coated on an ELISA plate. The unknown HNE protein samples or HNE-BSA standards are then added to the HNE conjugate preabsorbed ELISA plate. After a brief incubation, an anti-HNE polyclonal antibody is added, followed by an HRP conjugated secondary antibody. The content of HNE protein adducts in unknown samples is determined by comparison with a predetermined HNE-BSA standard curve.

Related Products

1. STA-305: OxiSelect™ Nitrotyrosine ELISA Kit
2. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit
3. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
4. STA-811: OxiSelect™ Methylglyoxal (MG) Competitive ELISA Kit
5. STA-813: OxiSelect™ N^ε-(carboxyethyl) lysine (CEL) Competitive ELISA Kit
6. STA-816: OxiSelect™ N^ε-(carboxymethyl) lysine (CML) Competitive ELISA Kit
7. STA-817: OxiSelect™ Advanced Glycation End Products (AGE) Competitive ELISA Kit
8. STA-832: OxiSelect™ MDA Adduct Competitive ELISA Kit

Kit Components

Box 1 (shipped at room temperature)

1. Protein Binding Strip Well Plate (Part No. 231001-T): One strip well microplate containing 32 wells (8 x 4).
2. Anti-HNE Antibody (1000X) (Part No. 283801-T): One 5 μ L vial of anti-HNE antibody.
3. Secondary Antibody, HRP Conjugate (1000X) (Part No. 231704): One 20 μ L vial.
4. Assay Diluent (Part No. 310804-T): One 20 mL bottle.
5. 10X Wash Buffer (Part No. 310806-T): One 30 mL bottle.
6. Substrate Solution (Part No. 310807-T): One 4 mL amber bottle.
7. Stop Solution (Part. No. 310808-T): One 4 mL bottle.

Box 2 (shipped on blue ice packs)

1. HNE-BSA Standard (Part No. 283803-T): One 100 μ L vial of 1 mg/mL HNE-BSA in PBS.
2. HNE Conjugate (Part No. 283802-T): One 20 μ L vial of HNE conjugate at 1.0 mg/mL in PBS.
3. 100X Conjugate Diluent (Part No. 281603-T): One 100 μ L vial.

Materials Not Supplied

1. Protein samples such as purified protein, plasma, serum, cell lysate
2. 1X PBS
3. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
4. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
5. Multichannel micropipette reservoir
6. Microplate reader capable of reading at 450 nm (620 nm as optional reference wave length)

Storage

Upon receipt, aliquot and store the Anti-HNE Antibody, HNE-BSA Standard, HNE Conjugate and 100X Conjugate Diluent at -20°C to avoid multiple freeze/thaw cycles. Store all other kit components at 4°C.

Preparation of Reagents

- HNE Conjugate Coated Plate:

Note: The HNE Conjugate coated wells are not stable and should be used within 24 hrs after coating. Only coat the number of wells to be used immediately.

1. Immediately before use, prepare 1X Conjugate Diluent by diluting the 100X Conjugate Diluent in 1X PBS. Example: Add 50 μ L to 4.95 mL of 1X PBS.

2. Immediately before use, prepare 10 µg/mL of HNE Conjugate by diluting the 1.0 mg/mL HNE Conjugate in 1X PBS. Example: Add 15 µL to 1.485 mL of 1X PBS.
 3. Mix the 10 µg/mL of HNE Conjugate and 1X Conjugate Diluent at 1:1 ratio and add 100 µL of the mixture to each well and incubate overnight at 4°C. Remove the HNE Conjugate coating solution and wash twice with 1X PBS. Blot plate on paper towels to remove excess fluid. Add 200 µL of Assay Diluent to each well and block for 1 hr at room temperature. Transfer the plate to 4°C and remove the Assay Diluent **immediately before use**.
- 1X Wash Buffer: Dilute the 10X Wash Buffer to 1X with deionized water. Stir to homogeneity.
 - Anti-HNE Antibody and Secondary Antibody: Immediately before use, dilute the Anti-HNE antibody 1:1000 and Secondary Antibody 1:1000 with Assay Diluent. Do not store diluted solutions.

Preparation of Standard Curve

Prepare a dilution series of HNE-BSA standards in the concentration range of 0 to 200 µg/mL by diluting the HNE-BSA Standard in Assay Diluent (Table 1).

Standard Tubes	1 mg/mL HNE-BSA Standard (µL)	Assay Diluent (µL)	HNE-BSA (µg/mL)
1	80	320	200
2	200 of Tube #1	200	100
3	200 of Tube #2	200	50
4	200 of Tube #3	200	25
5	200 of Tube #4	200	12.5
6	200 of Tube #5	200	6.25
7	200 of Tube #6	200	3.13
8	200 of Tube #7	200	1.56
9	0	200	0

Table 1. Preparation of HNE-BSA Standards

Assay Protocol

1. Prepare and mix all reagents thoroughly before use. Each HNE sample including unknown and standard should be assayed in duplicate.
2. Add 50 µL of unknown sample or HNE-BSA standard to the wells of the HNE Conjugate coated plate. If needed, unknown samples may be diluted in 1X PBS containing 0.1% BSA before adding. Incubate at room temperature for 10 minutes on an orbital shaker.
3. Add 50 µL of the diluted anti-HNE antibody to each well, incubate at room temperature for 1 hour on an orbital shaker.

4. Wash 3 times with 250 μ L of 1X Wash Buffer with thorough aspiration between each wash. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
5. Add 100 μ L of the diluted Secondary Antibody-HRP Conjugate to all wells and incubate for 1 hour at room temperature on an orbital shaker. Wash the strip wells 3 times according to step 4 above.
6. Warm Substrate Solution to room temperature. Add 100 μ L of Substrate Solution to each well. Incubate at room temperature for 2-20 minutes on an orbital shaker.
Note: Watch plate carefully; if color changes rapidly, the reaction may need to be stopped sooner to prevent saturation.
7. Stop the enzyme reaction by adding 100 μ L of Stop Solution to each well. Results should be read immediately (color will fade over time).
8. Read absorbance of each well on a microplate reader using 450 nm as the primary wave length.

Example of Results

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

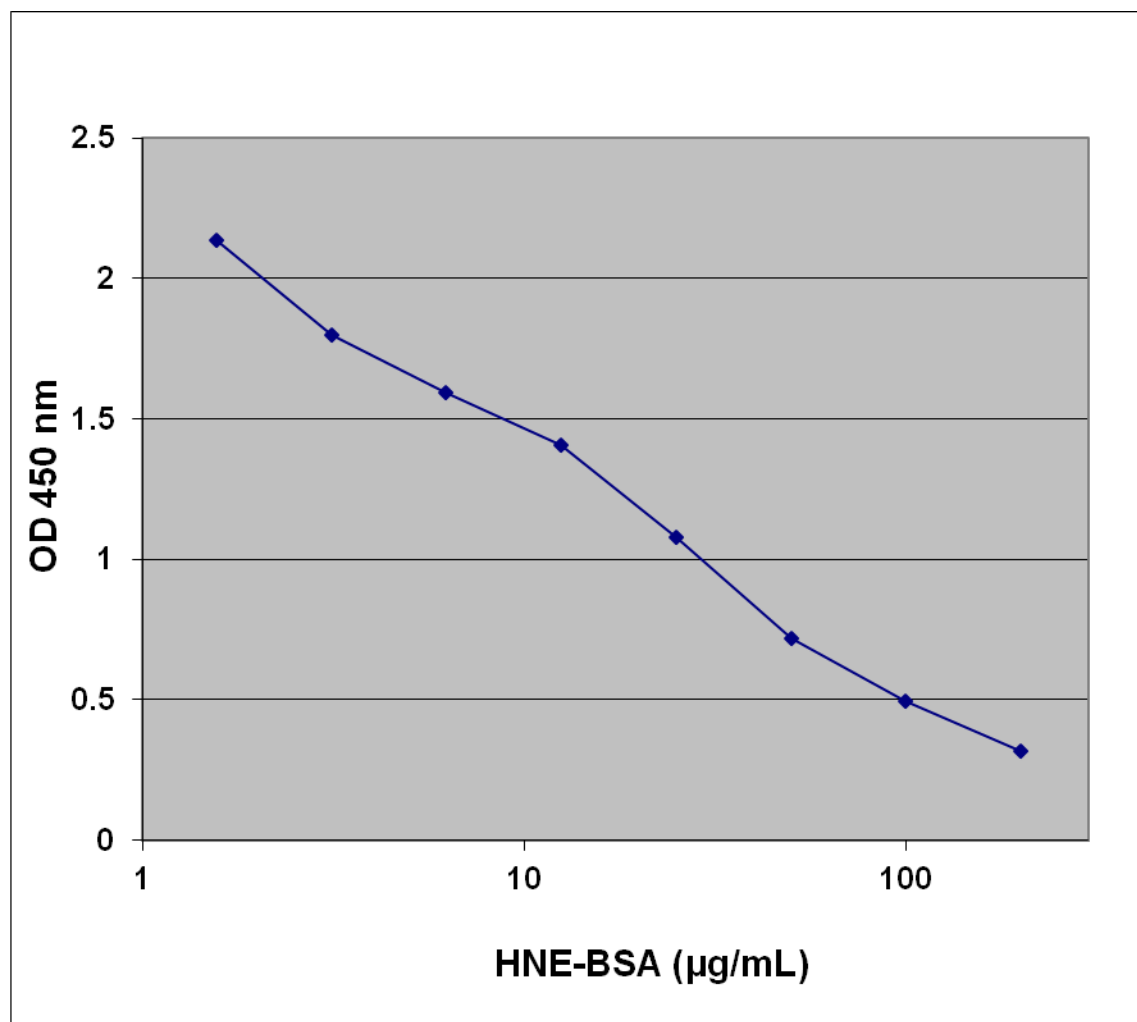


Figure 1: HNE-BSA Competitive ELISA Standard Curve.

References

1. Hoff HF, O'Neil J. (1993) *J Lipid Res.* 34: 1209-17.
2. Armstrong, D. and Browne, R. (1994). *Free Radicals in Diagnostic Medicine.* 366: 43-58.
3. Armstrong, D., et al. (1998). *Free Radicals and Antioxidant Protocols.* 108: 315-324.
4. Boyum, A. (1966). *J. of Clinical Investigation.* 21: Supplement 97.
5. Braun, D. and Fromherz, P. (1997). *Applied Physics A.*
6. Gidez, L., et al. (1982). *J. of Lipid Research.* 23: 1206-1223.
7. Lef'evre G., et al. (1998). *Annals de Biologie Clinique.* 56(3): 305-319.

8. Ohkawa, H., et al. (1979). *Anal. Biochem.* 95: 351-358.
9. Yagi, K. (1998). *Free Radicals and Antioxidant Protocols.* 108: 101-106.

Recent Product Citations

1. Kandimalla, R. et al. (2018). Hippocampal phosphorylated tau induced cognitive decline, dendritic spine loss and mitochondrial abnormalities in a mouse model of Alzheimer's disease. *Hum Mol Genet.* **27**(1):30-40. doi: 10.1093/hmg/ddx381.
2. Zhu, X. et al. (2017). Repeated restraint stress increases seizure susceptibility by activation of hippocampal endoplasmic reticulum stress. *Neurochemistry International.* **110**: 25-37.
3. Kołodziej, U. et al. (2017). Chronic high-protein diet induces oxidative stress and alters the salivary gland function in rats. *Arch Oral Biol.* **84**:6-12. doi: 10.1016/j.archoralbio.2017.09.006.
4. Bhowmick, S. and Drew, K.L. (2017). Arctic ground squirrel resist peroxynitrite-mediated cell death in response to oxygen glucose deprivation. *Free Radic Biol Med.* **113**:203-211. doi: 10.1016/j.freeradbiomed.2017.09.024.
5. Mazereeuw, G. et al. (2017). Baseline Oxidative Stress Is Associated with Memory Changes in Omega-3 Fatty Acid Treated Coronary Artery Disease Patients. *Cardiovasc Psychiatry Neurol.* 2017:3674371. doi: 10.1155/2017/3674371.
6. Tofiño-Vian, M. et al. (2017). Extracellular Vesicles from Adipose-Derived Mesenchymal Stem Cells Downregulate Senescence Features in Osteoarthritic Osteoblasts. *Oxid Med Cell Longev.* 2017:7197598. doi: 10.1155/2017/7197598.
7. Eraslan, G. et al. (2017). The effects of diosmin on aflatoxin-induced liver and kidney damage. *Environ Sci Pollut Res Int.* **24**(36):27931-27941. doi: 10.1007/s11356-017-0232-7.
8. Xu, W. et al (2017). MicroRNA-27b inhibition promotes Nrf2/ARE pathway activation and alleviates intracerebral hemorrhage-induced brain injury. *Oncotarget.* **8**(41):70669-70684. doi: 10.18632/oncotarget.19974. eCollection 2017 Sep 19.
9. Ren, Z. et al. (2017). Ethanol-induced Damage to the Developing Spinal Cord: The Involvement of CCR2 Signaling. *Biochim Biophys Acta.* pii: S0925-4439(17)30270-3. doi: 10.1016/j.bbadis.2017.07.035.
10. Meng, H. et al. (2017). Ameliorative Effect of Daidzein on Cisplatin-Induced Nephrotoxicity in Mice via Modulation of Inflammation, Oxidative Stress, and Cell Death. *Oxid Med Cell Longev.* 2017:3140680. doi: 10.1155/2017/3140680. Shen, K. et al. (2017). Baicalin ameliorates experimental liver cholestasis in mice by modulation of oxidative stress, inflammation, and NRF2 transcription factor. *Oxid. Med. Cell. Longev.* doi:10.1155/2017/6169128.
11. Shen, K. et al. (2017). Baicalin ameliorates experimental liver cholestasis in mice by modulation of oxidative stress, inflammation, and NRF2 transcription factor. *Oxid. Med. Cell. Longev.* doi:10.1155/2017/6169128.
12. Strickland, F.M. et al. (2017). Oxidative T cell modifications in lupus and Sjogren's syndrome. *Lupus* **2**(1):121.
13. Ouelaa, W. et al. (2017). Citrulline decreases hepatic endotoxin-induced injury in fructose-induced non-alcoholic liver disease: an ex vivo study in the isolated perfused rat liver. *Br. J. Nutr.* **22**:1-8.
14. Zhao, M. et al. Acrylamide-induced neurotoxicity in primary astrocytes and microglia: Roles of the Nrf2-ARE and NF-κB pathways. *Food Chem Toxicol.* **106**(Pt A):25-35.
15. Tahara, A. et al. (2017). Characterization and comparison of SGLT2 inhibitors: Part 3. Effects on diabetic complications in type 2 diabetic mice. *Eur. J. Pharmacol.* **809**:163-171.
16. Li, Q. et al (2017). Inhibition of neuronal ferroptosis protects hemorrhagic brain. *JCI Insight.* **2**:e90777. doi: 10.1172/jci.insight.90777.

17. Camaré, C. et al. (2017). 4-Hydroxynonenal Contributes to Angiogenesis through a Redox-Dependent Sphingolipid Pathway: Prevention by Hydralazine Derivatives. *Oxid Med Cell Longev.* 2017:9172741. doi: 10.1155/2017/9172741.
18. Pereira, I. et al. (2017). Transient Receptor Potential Ankyrin 1 Channel Expression on Peripheral Blood Leukocytes from Rheumatoid Arthritic Patients and Correlation with Pain and Disability. *Front Pharmacol.* **8**:53 doi: 10.3389/fphar.2017.00053.
19. Sussan, T.E. et al. (2017). Nrf2 regulates gene-environment interactions in an animal model of intrauterine inflammation: Implications for preterm birth and prematurity. *Sci Rep.* **7**:40194. doi: 10.1038/srep40194.
20. Verbeek, J. et al. (2017). Dietary intervention, but not losartan, completely reverses non-alcoholic steatohepatitis in obese and insulin resistant mice. *Lipids Health Dis.* **16**(1):46. doi: 10.1186/s12944-017-0432-7.
21. Young, B.E., et al. (2017). Markers of Oxidative Stress in Human Milk do not Differ by Maternal BMI But are Related to Infant Growth Trajectories. *Matern Child Health J.* doi:10.1007/s10995-016-2243-2
22. Velasco-Ortega E, et al. (2016). Dentistry and Diabetes: The Influence of Diabetes in Oral Diseases and Dental Treatments. *J Diabetes Res.* **2016**:6073190. doi: 10.1155/2016/6073190.
23. Newton, D.F. et al. (2016). Association of lipid peroxidation and brain-derived neurotrophic factor with executive function in adolescent bipolar disorder. *Psychopharmacology* doi:10.1007/s00213-016-4500-x.
24. Wang, X. et al. (2016). Fascin2 regulates cisplatin-induced apoptosis in NRK-52E cells. *Toxicol. Lett.* **266**:56-64.
25. Mukhopadhyay, P. et al. (2016). PARP inhibition protects against alcoholic and nonalcoholic steatohepatitis. *J. Hepatol.* doi:10.1016/j.jhep.2016.10.023.

Please see the complete list of product citations: <http://www.cellbiolabs.com/hne-adduct-competitive-elisa>.

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' 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

©2013-2018: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.