
Product Manual

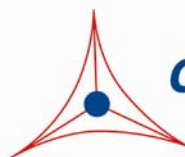
OxiSelect™ Protein Carbonyl Immunoblot Kit

Catalog Number

STA- 308

10 blots

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

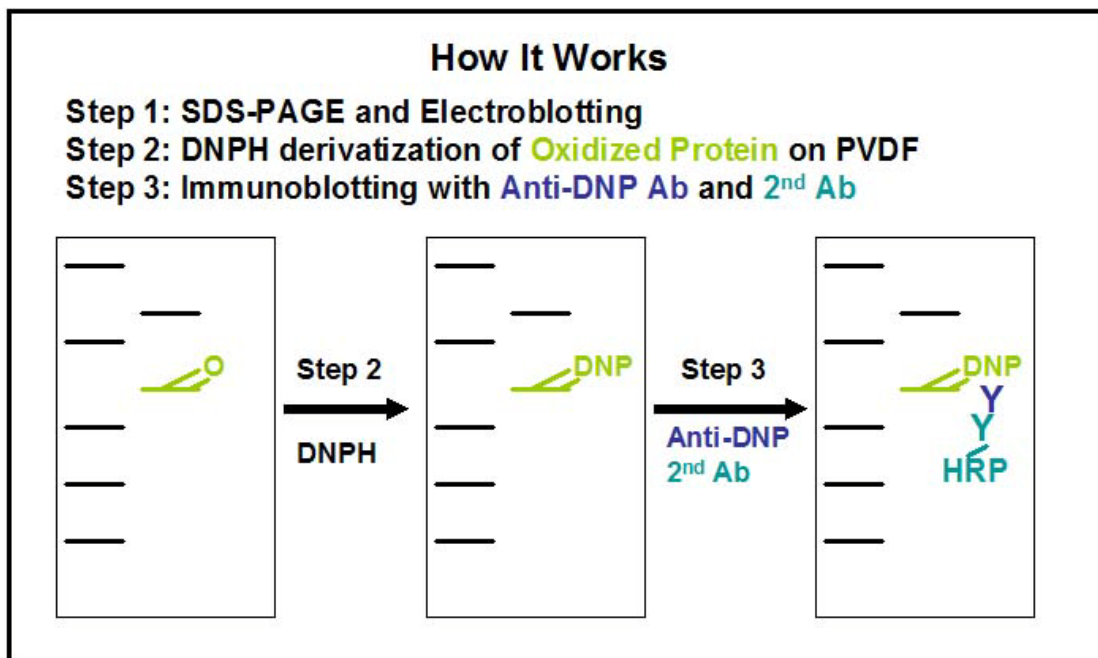
Protein oxidation is defined as the covalent modification of a protein induced either directly by reactive oxygen species or indirectly by reaction with secondary by-products of oxidative stress. Oxidative modification of proteins can be induced *in vitro* by a wide array of pro-oxidant agents and occurs *in vivo* during aging and in certain disease conditions.

There are numerous types of protein oxidative modification. The most common products of protein oxidation in biological samples are the protein carbonyl derivatives of Pro, Arg, Lys, and Thr. These derivatives are chemically stable and serve as markers of oxidative stress for most types of ROS.

Many of the current assays involve pre-derivatization of the carbonyl group with dinitrophenylhydrazine (DNPH) prior to electrophoresis, followed by immunoblotting with an anti-DNP antibody. Unfortunately, this pre-derivatization alters the electrophoretic (and electrofocusing) properties of proteins. Consequently, it is not possible to directly compare the patterns from "oxidized" fingerprints with those from "non-oxidized" protein fingerprints. Cell Biolabs' Protein Carbonyl Immunoblot Kit has the ability to conduct all derivatization and staining **after** electrophoresis and transblotting. This allows one to directly compare oxidized vs. non-oxidized protein fingerprints.

The OxiSelect™ Protein Carbonyl Immunoblot Kit offers a simple and complete system for the detection of protein oxidation. This kit also includes a Protein Oxidation Immunoblot Control as positive control. Each kit provides sufficient quantities to perform at least 10 blots (7.5 cm X 8.5 cm).

Assay Principle



Related Products

1. STA-303: OxiSelect™ Nitrotyrosine Protein Immunoblot Kit
2. STA-304: Protein Tyrosine Nitration Control (Nitrotyrosine-BSA)
3. STA-305: OxiSelect™ Nitrotyrosine ELISA Kit
4. STA-309: Oxidized Protein Immunoblot Control (Carbonyl-BSA)
5. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit
6. STA-315: OxiSelect™ Protein Carbonyl Spectrophotometric Assay
7. STA-318: OxiSelect™ AOPP Assay Kit
8. STA-319: AOPP-Human Serum Albumin (AOPP-HSA)
9. STA-816: OxiSelect™ N-epsilon-(Carboxymethyl) Lysine (CML) Competitive ELISA Kit
10. STA-817: OxiSelect™ Advanced Glycation End Products (AGE) Competitive ELISA

Kit Components

1. Rabbit Anti-DNP Antibody (Part No. 230801): One 100 µL tube
2. Secondary Antibody, HRP-conjugate (Part No. 230805): One 100 µL tube
3. Protein Oxidation Immunoblot Control (Part No. 230803): One 100 µL tube (provided ready-to-use oxidized BSA in 1X reducing SDS-PAGE Sample Buffer, pre-boiled)
4. 10X DNPH Solution (Part No. 230804): One 20 mL amber bottle

Materials Not Supplied

1. Protein MW standards
2. Reducing SDS-PAGE Sample Buffer
3. Polyacrylamide gels such as precast gels available from Invitrogen or BioRad
4. Electrophoresis Buffers
5. Electrophoresis and Western Blot Transfer Systems
6. Immunoblotting Buffers such as TBST (20 mM Tris-HCl, pH 7.4, 0.15 M NaCl, 0.05% Tween-20)
7. PVDF or Nitrocellulose Membrane (PVDF is recommended)
8. Methanol
9. 2N HCl
10. Non-fat Dry Milk
11. ECL Reagents

Storage

Upon receipt, store the 10X DNPH Solution at 4°C. Aliquot and store all other components at -20°C to avoid multiple freeze/thaw cycles.

Assay Protocol

I. Electrophoresis and Transblotting

1. Prepare samples for electrophoresis with reducing SDS Sample Buffer.
2. Load 20 µL of Protein Oxidation Immunoblot Control (provided ready-to-use, pre-boiled) or sample to wells of a polyacrylamide gel. Also, it's recommended to include a pre-stained MW standard (as indicator of a successful transfer in step 3). Run the gel as per the manufacturer's instructions.
3. Transfer the gel proteins to a PVDF membrane as per the manufacturer's instructions.

Note: We recommend using PVDF membrane instead of Nitrocellulose due to its low background signal after derivatization, resulting in stronger chemiluminescent signal.

II. Derivatization (all steps are at room temperature, with shaking)

1. Following the electroblotting step, immerse the PVDF membrane in 100% Methanol for 15 seconds, and then allow it to dry at room temperature for 5 minutes.

Note: If Nitrocellulose is used instead of PVDF, this step should be skipped.

2. Equilibrate the membrane in TBS containing 20% Methanol for 5 minutes.
3. Wash the membrane in 2N HCl for 5 minutes.
4. Prepare sufficient amount of 1X DNPH solution by diluting the 10X DNPH Solution in 2N HCl. Incubate the membrane with 1X DNPH solution for exactly 5 minutes.

Note: 1X DNPH Solution is stable for one week when stored in the dark at 4°C. Do not freeze.

5. Wash the membrane three times in 2N HCl, 5 minutes each time.
6. Next, wash the membrane five times in 100% methanol (PVDF) or 50% methanol (Nitrocellulose), 5 minutes each time.

III. Immunoblotting

1. Block the DNPH-treated membrane with 5% non-fat dry milk in TBST for 1 hr at room temperature with constant agitation.
2. Wash the blocked membrane three times with TBST, 5 minutes each time.
3. Incubate the membrane with Rabbit Anti-DNP Antibody, freshly diluted 1:1000 in 5% non-fat dry milk/TBST, for 1-2 hr at room temperature with constant agitation.
4. Wash the blotted membrane three times with TBST, 5 minutes each time.
5. Incubate the membrane with Secondary Antibody, HRP-conjugate, freshly diluted 1:1000 in 5% non-fat dry milk/TBST, for 1 hr at room temperature with constant agitation.

6. Wash the blotted membrane five times with TBST, 5 minutes each time.
7. Use the detection method of your choice. We recommend enhanced chemiluminescence reagents from Pierce.

Example of Results

The following figure demonstrates typical blot results of oxidized BSA after DNPH derivation. One should use the data below for reference only. This data should not be used to interpret actual results.



Figure 1: Immunoblotting of Oxidized BSA. Carbonyl-BSA, Oxidation Immunoblot Control, was first electroblotted onto nitrocellulose membrane. Following the electroblotting procedure, the membrane was incubated with (right strip) or without (left strip) DNPH solution. The derivatized Carbonyl-BSA is detected by immunoblotting with anti-DNP antibody as described in the Assay Protocol.

References

1. Cadenas, E., Boveris, A., Ragan, CI., and Stoppani, AO. (1977). *Archives of Biochemistry & Biophysics* **180**:248–257.
2. Wakeyama, H., Takeshige, K., Takayanagi, R., and Minakami, S. (1982). *Biochem J.* **205**:593–601.
3. Talent, JM., Kong, Y., and Gracy, RW. (1998). *Anal. Biochem.* **263**:31–38.

Recent Product Citations

1. Li, Y. et al. (2017). Changes in the mitochondrial protein profile due to ROS eruption during ageing of elm (*Ulmus pumila* L.) seeds. *Plant Physiol Biochem.* **114**:72-87. doi: 10.1016/j.plaphy.2017.02.023.

2. Konopka, A.R. et al. (2017). Influence of Nrf2 activators on subcellular skeletal muscle protein and DNA synthesis rates after 6 weeks of milk protein feeding in older adults. *Geroscience*. doi: 10.1007/s11357-017-9968-8.
3. Kazi, R.S. et al. (2017). Glycation inhibitors extend yeast chronological lifespan by reducing advanced glycation end products and by back regulation of proteins involved in mitochondrial respiration. *J Proteomics*. **156**:104-112. doi: 10.1016/j.jprot.2017.01.015.
4. Galmes-Pascual, B.M. et al. (2016). 17 β -estradiol improves hepatic mitochondrial biogenesis and function through PGC1B. *J. Endocrinol.* **232**:297-308.
5. Su, L. et al. (2016). Reactive oxygen species induced by cold stratification promote germination of *Hedysarum scoparium* seeds. *Plant Physiol. Biochem.* **109**:406-415.
6. Ferrer, M.D. et al. (2016). Haem biosynthesis and antioxidant enzymes in circulating cells of acute intermittent porphyria patients. *PLoS One* doi:10.1371/journal.pone.0164857.
7. Capo, X. et al. (2016). Effects of dietary almond-and olive oil-based docosahexaenoic- and vitamin E-enriched beverage supplementation on athletic performance and oxidative stress. *Food Funct.* **7**:4920-4934.
8. Chaves, R. S. et al. (2016). Presence of insoluble Tau following rotenone exposure ameliorates basic pathways associated with neurodegeneration. *IBRO Rep.* doi:10.1016/j.ibror.2016.09.001.
9. Almeida, M. F. et al.. Aged Lewis rats exposed to low and moderate doses of rotenone are a good model for studying the process of protein aggregation and its effects upon central nervous system cell physiology. *Arq Neuropsiquiatr.* **74**:737-744.
10. Ong, L. K. et al. (2016). Reconsidering the role of glial cells in chronic stress-induced dopaminergic neurons loss within the substantia nigra? Friend or foe? *Brain Behav Immun.* doi:10.1016/j.bbi.2016.10.001.
11. Busquets-Cortés, C. et al. (2016). Training enhances immune cells mitochondrial biosynthesis, fission, fusion, and their antioxidant capabilities synergistically with dietary docosahexaenoic supplementation. *Oxid Med Cell Longev.* doi:10.1155/2016/8950384.
12. Hudgens, J. L. et al. (2016). Platelet-rich plasma activates proinflammatory signaling pathways and induces oxidative stress in tendon fibroblasts. *Am J Sports Med.* doi:10.1177/0363546516637176.
13. Laitano, O. et al. (2016). Pharmacological targeting of mitochondrial reactive oxygen species counteracts diaphragm weakness in chronic heart failure. *J Appl Physiol.* doi:10.1152/jappphysiol.00822.2015.
14. Tanase, M. et al. (2016). Role of carbonyl modifications on aging-associated protein aggregation. *Sci Rep.* doi:10.1038/srep19311.
15. ElHajj, Z. et al. (2016). Effects of postmortem delays on protein composition and oxidation. *Brain Res Bull.* doi:10.1016/j.brainresbull.2016.01.005.
16. Sbert-Roig, M. et al. (2015). GPER mediates the effects of 17 β -estradiol in cardiac mitochondrial biogenesis and function. *Mol Cell Endocrinol.* doi:10.1016/j.mce.2015.11.027.
17. Zhou, J. et al. (2015). Correlations between photodegradation of bisretinoid constituents of retina and dicarbonyl adduct deposition. *J Biol Chem.* **290**:27215-27227.
18. Martorell, M. et al. (2015). Docosahexaenoic acid supplementation promotes erythrocyte antioxidant defense and reduces protein nitrosative damage in male athletes. *Lipids.* **50**:131-148.
19. Capó, X. et al. (2015). Diet supplementation with DHA-enriched food in football players during training season enhances the mitochondrial antioxidant capabilities in blood mononuclear cells. *Eur J Nutr.* **54**:35-49.
20. Cui, J. et al. (2015). Rapamycin protects against gentamicin-induced acute kidney injury via autophagy in mini-pig models. *Sci Rep.* doi:10.1038/srep11256.

21. Konopka, A. R. et al. (2015). Defects in mitochondrial efficiency and H₂O₂ emissions in obese women are restored to a lean phenotype with aerobic exercise training. *Diabetes*. **64**:2104-2115.
22. Pons, D. G. et al. (2015). UCP2 inhibition sensitizes breast cancer cells to therapeutic agents by increasing oxidative stress. *Free Radic Biol Med*. doi:10.1016/j.freeradbiomed.2015.04.032.
23. Nadal-Serrano, M. et al. (2015). Chronic-leptin attenuates cisplatin cytotoxicity in MCF-7 breast cancer cell line. *Cell Physiol Biochem*. **36**:221-232.
24. Blanquer-Rosselló, M. D. et al. (2015). Leptin modulates mitochondrial function, dynamics and biogenesis in MCF-7 cells. *J Cell Biochem*. doi: 10.1002/jcb.25158. Cao, W. et al. (2015). Geranylgeranylacetone ameliorates lung ischemia/reperfusion injury by HSP70 and thioredoxin redox system: NF-κB pathway involved. *Pulm Pharmacol Ther*. doi: 10.1016/j.pupt.2015.02.009.
25. Konopka, A. R. et al. (2015). Defects in mitochondrial efficiency and H₂O₂ emissions in obese women are restored to a lean phenotype with aerobic exercise training. *Diabetes* doi: 10.2337/db14-1701.
26. Martorell, M. et al. (2014). Docosahexaenoic acid supplementation promotes erythrocyte antioxidant defense and reduces protein nitrosative damage in male athletes. *Lipids*. **50**:131-148.
27. Wang, Z. H. et al. (2014). Concentration-dependent wrestling between detrimental and protective effects of H₂O₂ during myocardial ischemia/reperfusion. *Cell Death Dis*. **5**:e1297.
28. Chen, C. et al. (2014). C75, An inhibitor of fatty acid synthase, suppresses the mitochondrial fatty acid synthesis pathway and impairs mitochondrial function. *J Biol Chem*. **289**:17184-17194.
29. Hwee, D. T. et al. (2014). Maintenance of muscle mass and load-induced growth in Muscle RING Finger 1 null mice with age. *Aging Cell*. **13**:92-101.
30. Galay, R. L. et al. (2014). Two kinds of ferritin protect ixodid ticks from iron overload and consequent oxidative stress. *PLoS One*. **9**:e90661.

Please see the complete list of product citations: <http://www.cellbiolabs.com/protein-carbonyl-immunoblot-kit>

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

©2004-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.