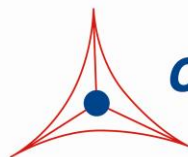

Product Manual

OxiSelect™ Nitrotyrosine ELISA Kit

Catalog Number

STA-305	96 assays
STA-305-5	5 x 96 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

The modification of tyrosine residues in proteins to 3-nitrotyrosine by peroxynitrite (Figure 1) or other potential nitrating agents has been detected in biological systems that are subject to oxidative stress. Detection of nitrotyrosine-containing proteins has been reported in many human and animal diseases or cellular models of disease. While all tyrosine residues in proteins may theoretically be targets for nitration, presumably the efficiency of tyrosine nitration is dependent on various biological conditions such as the local production and concentration of the reactive species, the existence and availability of antioxidants and scavengers, the accumulation of inflammatory cell and the presence of pro-inflammatory cytokines, as well as the proximity and compartmentation of these components.

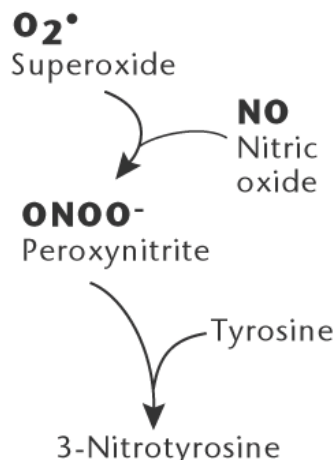


Figure 1. 3-Nitrotyrosine Formation

The OxiSelect™ Nitrotyrosine ELISA Kit is a competitive enzyme immunoassay developed for rapid detection and quantitation of 3-nitrotyrosine in protein sample. The quantity of 3-nitrotyrosine in protein sample is determined by comparing its absorbance with that of a known nitrated BSA standard curve. The kit has a nitrotyrosine detection sensitivity range of 20 nM to 8.0 μ M. Each kit provides sufficient reagents to perform up to 96 assays, including standard curve and unknown protein samples.

Assay Principle

The nitrotyrosine quantitation kit is a competitive ELISA. The unknown protein nitrotyrosine sample or nitrated BSA standards are first added to a nitrated BSA preabsorbed EIA plate. After a brief incubation, an anti-nitrotyrosine antibody is added, followed by an HRP conjugated secondary antibody. The protein nitrotyrosine content in unknown sample is determined by comparing with a standard curve that is prepared from predetermined nitrated BSA standards.

Related Products

1. STA-304: Protein Tyrosine Nitration Control (Nitrotyrosine-BSA)
2. STA-308: OxiSelect™ Protein Carbonyl Immunoblot Kit
3. STA-309: Oxidized Protein Immunoblot Control (Carbonyl-BSA)
4. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit

5. STA-318: OxiSelect™ AOPP Assay Kit
6. STA-319: AOPP-Human Serum Albumin (AOPP-HSA)
7. STA-816: OxiSelect™ N-epsilon-(Carboxymethyl) Lysine (CML) Competitive ELISA Kit
8. STA-817: OxiSelect™ Advanced Glycation End Products (AGE) Competitive ELISA Kit

Kit Components

1. Nitrotyrosine Coated EIA Plate (Part No. 230501): One strip well 96-well plate.
2. Anti-Nitrotyrosine Antibody (Part No. 230502): One 20 µL vial of anti-nitrotyrosine Rabbit IgG.
3. Secondary Antibody, HRP Conjugate (Part No. 231009): One 20 µL 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.
8. Nitrated BSA Standard (Part No. 230503): One 500 µL vial of 1 mg/mL Nitrated BSA in PBS with a nitrotyrosine content of 40 µM (2.7 mole of nitrotyrosine per mole of BSA). The protein nitrotyrosine level is predetermined by a spectrophotometric method as described by Ischiropoulos et al (See Ref. 3).

Materials Not Supplied

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

Storage

Upon receipt, aliquot and store the Nitrated BSA Standard at -20°C to avoid multiple freeze/thaw cycles. Store all other kit components at 4°C.

Preparation of Reagents

- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.
- Anti-Nitrotyrosine Antibody and Secondary Antibody: Immediately before use dilute the Anti-Nitrotyrosine 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 Nitrated BSA Standards in the nitrotyrosine concentration range of 0 nM – 8000 nM by diluting the Nitrated BSA stock solution in Assay Diluent (Table 1).

Standard Tubes	Nitrated BSA Standard (µL)	Assay Diluent (µL)	Nitrated BSA (µg/mL)	Nitrotyrosine (nM)
1	60	240	200	8000
2	100 of Tube #1	300	50	2000
3	100 of Tube #2	300	12.5	500
4	100 of Tube #3	300	3.125	125
5	100 of Tube #4	300	0.78	31.25
6	100 of Tube #5	300	0.195	7.81
7	100 of Tube #6	300	0.049	1.95
8	0	300	0	0

Table 1. Preparation of Nitrated BSA Standards

Assay Protocol

1. Prepare and mix all reagents thoroughly before use. Each protein sample including nitrated BSA and blank should be assayed in duplicate.
2. Add 50 µL of unknown protein sample or nitrated BSA standard to the wells of the EIA plate. Incubate at room temperature for 10 minutes on an orbital shaker.
3. Add 50 µL of the diluted anti-nitrotyrosine antibody to each well, incubate at room temperature for 1 hour on an orbital shaker.
4. Wash microwell strips 3 times with 250 µL 1X Wash Buffer per well 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 µL of the diluted Secondary Antibody-Enzyme Conjugate to all wells.
6. Incubate at room temperature for 1 hour on an orbital shaker.
7. Wash microwell strips 3 times according to step 4 above. Proceed immediately to the next step.
8. Warm Substrate Solution to room temperature. 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 Nitrotyrosine ELISA results. One should use the data below for reference only. This data should not be used to interpret actual results.

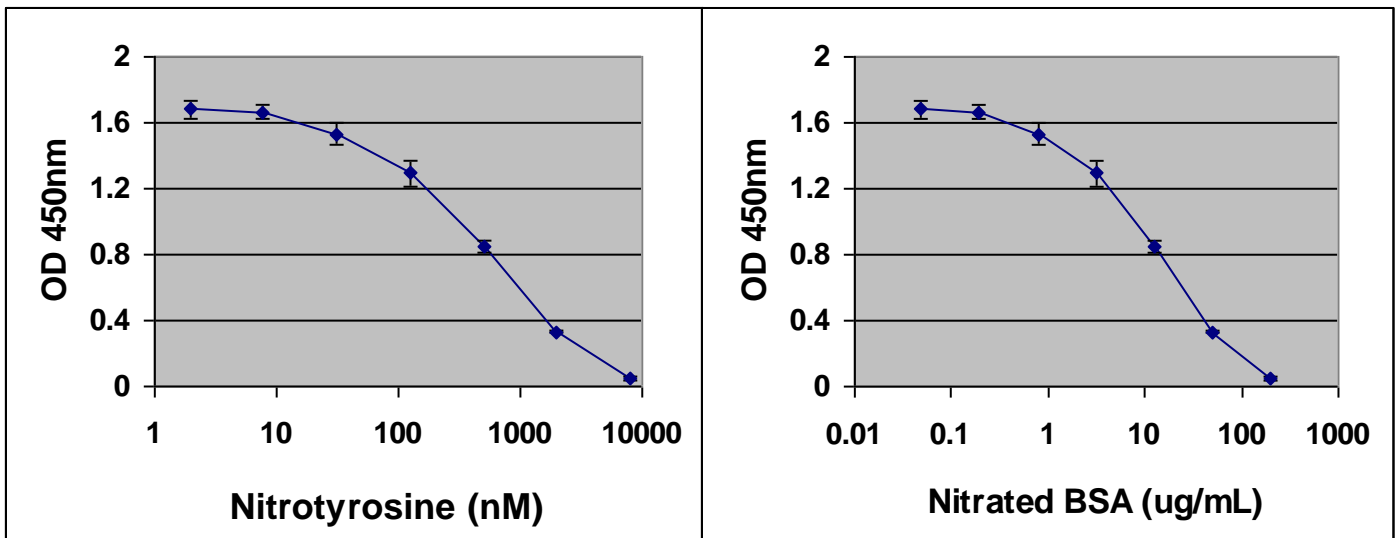


Figure 2: Nitrotyrosine ELISA Standard Curve.

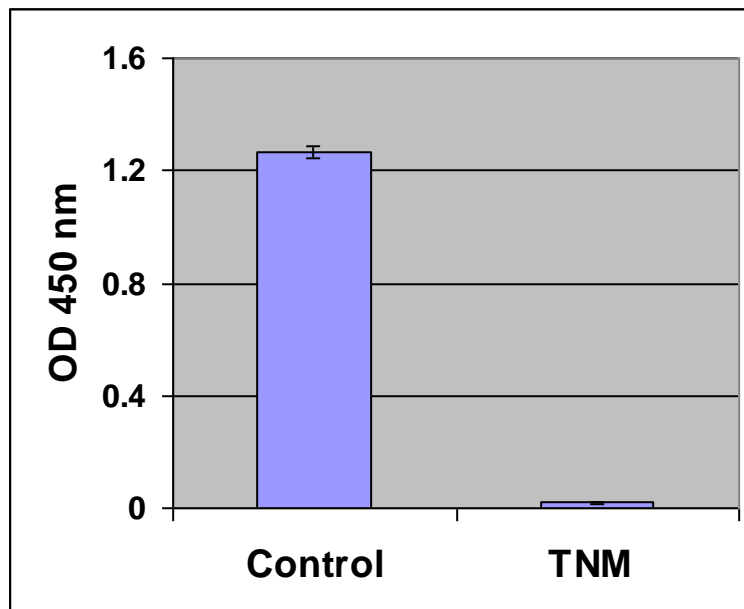


Figure 3: Protein Nitration by tetranitromethane. STO (MEF) cells were lysed in 25mM HEPES, pH 7.5, 150 mM NaCl, 1% NP-40, 10 mM MgCl₂, 1 mM EDTA, 2% Glycerol. Cell Lysate was nitrated with tetranitromethane (TNM). The protein 3-nitrotyrosone levels were determined as described in the assay instructions.

References

1. Gow A. J., Farkouh C. R., Munson D. A., Posencheg M. A and Ischiropoulos H. (2004) Am J Physiol Lung Cell Mol Physiol. **287**, L262-8.
2. Bian, K., Ke, Y., Kamisaki, Y. and Murad, F. (2006) J Pharmacol Sci Vol. **101**, 271-279.

3. Ischiropoulos, H., Zhu, L., Chen, J., Tsai, M., Martin, J. C., Smith, C. D. and Beckman, J. S. (1992) *Archiv. Biochem. Biophys.* **298**, 431-437.

Recent Product Citations

1. Maciejczyk, M. et al. (2022). α -Lipoic Acid Strengthens the Antioxidant Barrier and Reduces Oxidative, Nitrosative, and Glycative Damage, as well as Inhibits Inflammation and Apoptosis in the Hypothalamus but Not in the Cerebral Cortex of Insulin-Resistant Rats. *Oxid Med Cell Longev.* doi: 10.1155/2022/7450514.
2. Rosa, R.H. et al. (2022). Intravitreal Administration of Stanniocalcin-1 Rescues Photoreceptor Degeneration with Reduced Oxidative Stress and Inflammation in a Porcine Model of Retinitis Pigmentosa. *Am J Ophthalmol.* doi: 10.1016/j.ajo.2022.03.014.
3. Rodríguez-Pérez, M.D. et al. (2022). Neuroprotective Effect of 3',4'-Dihydroxyphenylglycol in Type-1-like Diabetic Rats-Influence of the Hydroxytyrosol/3',4'-dihydroxyphenylglycol Ratio. *Nutrients.* **14**(6):1146. doi: 10.3390/nu14061146.
4. Kosutova, P. et al. (2021). Nitric-Oxide-Releasing Dexamethasone Derivative NCX-1005 Improves Lung Function and Attenuates Inflammation in Experimental Lavage-Induced ARDS. *Pharmaceutics.* **13**(12):2092. doi: 10.3390/pharmaceutics13122092.
5. Williamson-Reisdorph, C.M. et al. (2021). Cardiovascular and Blood Oxidative Stress Responses to Exercise and Acute Woodsmoke Exposure in Recreationally Active Individuals. *Wilderness Environ Med.* doi: 10.1016/j.wem.2021.10.002.
6. Murata, T. et al. (2021). Chemical inducer of regucalcin attenuates lipopolysaccharide-induced inflammatory responses in pancreatic MIN6 β -cells and RAW264.7 macrophages. *FEBS Open Bio.* doi: 10.1002/2211-5463.13321.
7. Całyniuk, Z. et al. (2021). Selected metabolic, epigenetic, nitration and redox parameters in turkeys fed diets with different levels of arginine and methionine. *Ann. Anim. Sci.* doi: 10.2478/aoas-2021-0069.
8. Ognik, K. et al. (2021). The immune status, oxidative and epigenetic changes in tissues of turkeys fed diets with different ratios of arginine and lysine. *Sci Rep.* **11**(1):15975. doi: 10.1038/s41598-021-95529-y.
9. Williamson-Reisdorph, C.M. et al. (2021). Blood oxidative stress and post-exercise recovery are unaffected by hypobaric and hypoxic environments. *J Sports Sci.* doi: 10.1080/02640414.2021.1872960.
10. Baker, B.C. et al. (2021). Hypoxia and oxidative stress induce sterile placental inflammation in vitro. *Sci Rep.* **11**(1):7281. doi: 10.1038/s41598-021-86268-1.
11. Šutulović, N. et al. (2021). Experimental Chronic Prostatitis/Chronic Pelvic Pain Syndrome Increases Anxiety-Like Behavior: The Role of Brain Oxidative Stress, Serum Corticosterone, and Hippocampal Parvalbumin-Positive Interneurons. *Oxid Med Cell Longev.* doi: 10.1155/2021/6687493.
12. Jankowski, J. et al. (2021). The effect of different dietary ratios of lysine, arginine and methionine on protein nitration and oxidation reactions in turkey tissues and DNA. *Animal.* doi: 10.1016/j.animal.2021.100183.
13. Dietrich-Muszalska, A. et al. (2021). Comparative Study of the Effects of Atypical Antipsychotic Drugs on Plasma and Urine Biomarkers of Oxidative Stress in Schizophrenic Patients. *Neuropsychiatr Dis Treat.* **17**:555-565. doi: 10.2147/NDT.S283395.

14. Zhou, Q. et al. (2021). Physiological Responses of *Microcystis aeruginosa* to Extracellular Degradative Enzymes and Algicidal Substance from Heterotrophic Bacteria. *Pol. J. Environ. Stud.* **30**(3): 1-9. doi: 10.15244/pjoes/127867.
15. Ohira, H. et al. (2021). Alteration of oxidative-stress and related marker levels in mouse colonic tissues and fecal microbiota structures with chronic ethanol administration: Implications for the pathogenesis of ethanol-related colorectal cancer. *PLoS One.* **16**(2):e0246580. doi: 10.1371/journal.pone.0246580.
16. Uzunköprü, C. et al. (2021). Retinal Nerve Fiber Layer Thickness Correlates with Serum and Cerebrospinal Fluid Neurofilament Levels and is Associated with Current Disability in Multiple Sclerosis. *Arch Neuropsychiatry.* doi: 10.29399/npa.27355.
17. Gerszi, D. et al. (2021). Evaluation of oxidative/nitrative stress and uterine artery pulsatility index in early pregnancy. *Physiol Int.* doi: 10.1556/2060.2020.00041.
18. Wong, T.H.T. et al. (2020). Consuming decaffeinated coffee with milk and sugar added before a high-glycaemic index meal improves postprandial glycaemic and insulinaemic responses in healthy adults. *Br J Nutr.* **124**(8):785-796. doi: 10.1017/S0007114520001750.
19. Gil, A. et al. (2020). Neuronal Metabolism and Neuroprotection: Neuroprotective Effect of Fingolimod on Menadione-Induced Mitochondrial Damage. *Cells.* **10**(1):E34. doi: 10.3390/cells10010034.
20. Miller, V.J. et al. (2020). A Ketogenic Diet Combined With Exercise Alters Mitochondrial Function In Human Skeletal Muscle While Improving Metabolic Health. *Am J Physiol Endocrinol Metab.* doi: 10.1152/ajpendo.00305.2020.
21. Sprick, J.D. et al. (2020). Augmented exercise pressor response during maximal treadmill exercise is not related to systemic inflammation in stroke survivors. *Top Stroke Rehabil.* doi: 10.1080/10749357.2020.1806436.
22. Hamidi, V. et al. (2020). Acute Exenatide Therapy Attenuates Postprandial Vasodilation in Humans with Prediabetes: A Randomized Controlled Trial. *Metab Syndr Relat Disord.* doi: 10.1089/met.2019.0102.
23. Valvassori, S.S. et al. (2020). Protein kinase C isoforms as a target for manic-like behaviors and oxidative stress in a dopaminergic animal model of mania. *Prog Neuropsychopharmacol Biol Psychiatry.* doi: 10.1016/j.pnpbp.2020.109940.
24. Vujacic-Mirski, K. et al. (2020). Development of an Analytical Assay for Electrochemical Detection and Quantification of Protein-Bound 3-Nitrotyrosine in Biological Samples and Comparison with Classical, Antibody-Based Methods. *Antioxidants (Basel).* **9**(5). pii: E388. doi: 10.3390/antiox9050388.
25. Chira, A. et al. (2020). Monocyte chemotactic protein-1 and nitrotyrosine in irritable bowel syndrome. *Exp Ther Med.* doi: 10.3892/etm.2020.8665.
26. Casós, K. et al. (2020). Determination of Redox Status in Serum. *Methods Mol Biol.* **2110**:115-128. doi: 10.1007/978-1-0716-0255-3_8.
27. Ruottinen, M. et al. (2020). The Rectus Sheath Block (RSB) Analgesia Following Laparotomy Could Affect Malondialdehyde (MDA) Concentrations in Benign Disease and Cancer. *Anticancer Res.* **40**(1):253-259. doi: 10.21873/anticancer.13947.
28. Menegas, S. et al. (2020). Efficacy of folic acid as an adjunct to lithium therapy on manic-like behaviors, oxidative stress and inflammatory parameters in an animal model of mania. *Metab Brain Dis.* doi: 10.1007/s11011-019-00503-3.

29. Mikolka, P. et al. (2019). Effect of different dosages of dexamethasone therapy on lung function and inflammation in an early phase of acute respiratory distress syndrome model. *Physiol Res.* **68**(Supplementum 3):S253-S263. doi: 10.33549/physiolres.934364.
30. Balasubramanian, P. et al. (2019). Obesity-Induced Sympathoexcitation is Associated with Nrf2 Dysfunction in the Rostral Ventrolateral Medulla. *Am J Physiol Regul Integr Comp Physiol.* doi: 10.1152/ajpregu.00206.2019.

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

©2013-2022: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.