
Product Manual

OxiSelect™ Nitrotyrosine ELISA Kit, Trial Size

Catalog Number

STA-305-T

32 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.

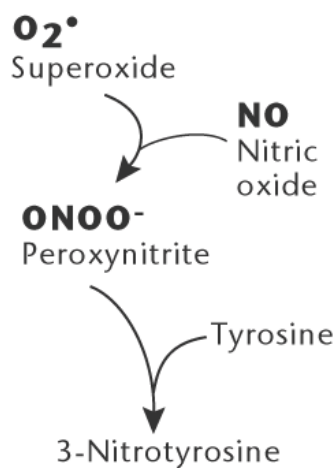


Figure 1. 3-Nitrotyrosine Formation

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 μM. Each Trial Size Nitrotyrosine ELISA Kit provides sufficient reagents to perform up to 32 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-303: OxiSelect™ Nitrotyrosine Protein Immunoblot Kit
2. STA-304: Protein Tyrosine Nitration Control (Nitrotyrosine-BSA)

3. STA-308: OxiSelect™ Protein Carbonyl Immunoblot 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

Kit Components

1. Nitrotyrosine Coated EIA Plate (Part No. 230501-T): One strip-well microplate containing 32 wells (8 x 4).
2. Anti-Nitrotyrosine Antibody (Part No. 230502-T): One 5 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-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.
8. Nitrated BSA Standard (Part No. 230503-T): One 200 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 microplate. 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.

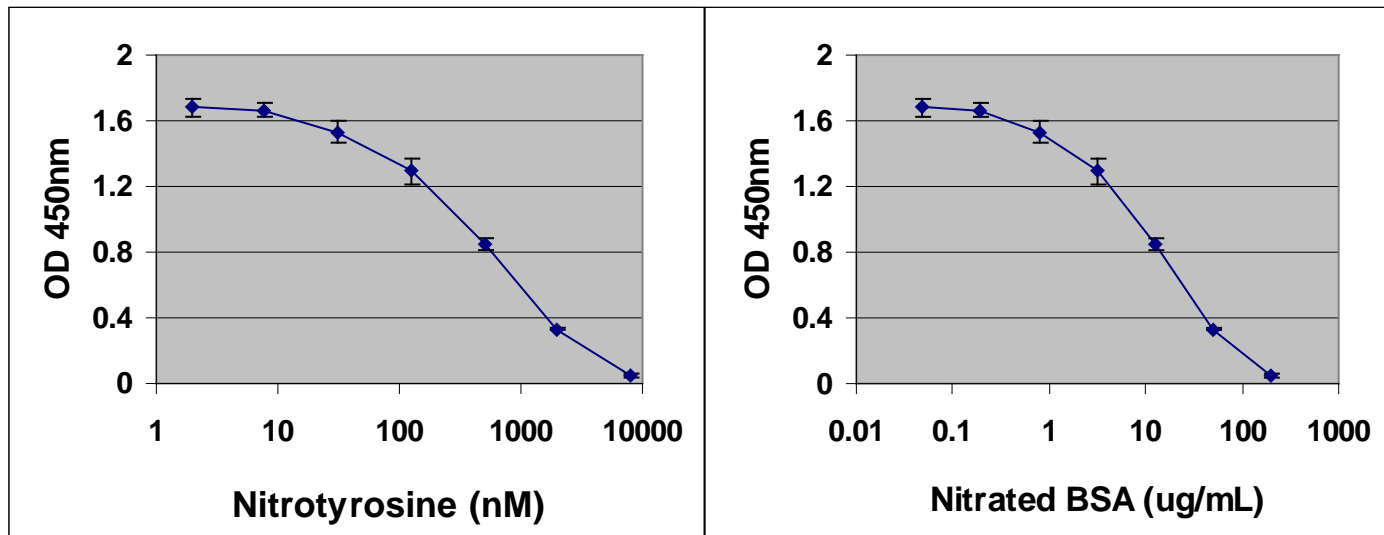


Figure 2: Nitrotyrosine ELISA Standard Curve.

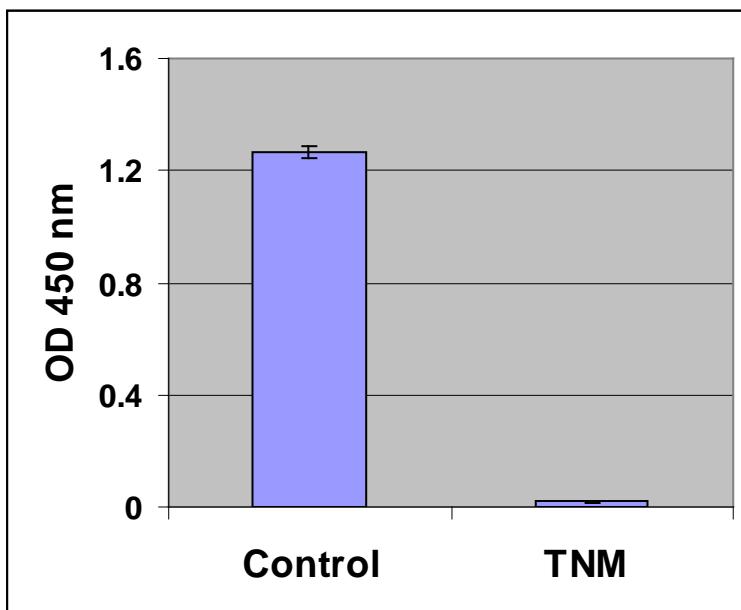


Figure 3: Protein Nitration by tetranitromethane. STO (MEF) cells were lysed in 25mM HEPES, pH 7.5, 150 mM NaCl, 1% NP-40, 10 mM MgCl₂, 1 mM EDTA, 2% Glycerol. Cell Lysate was nitrated with tetranitromethane (TNM). The protein 3-nitrotyrsonne levels were determined as described in the assay instructions.

References

1. Montez, P. et al. (2012). Angiotensin Receptor Blockade Recovers Hepatic UCP2 Expression and Aconitase and SDH Activities and Ameliorates Hepatic Oxidative Damage in Insulin Resistant Rats. *Endocrinology*. **153**:5845-5856
2. 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.
3. Bian, K., Ke, Y., Kamisaki, Y. and Murad, F. (2006) J Pharmacol Sci Vol. **101**, 271-279.
4. 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. Wan P, et al. (2017). Trimetazidine protects retinal ganglion cells from acute glaucoma via the Nrf2/Ho-1 pathway. *Clin Sci (Lond)*. **131**(18):2363-2375. doi: 10.1042/CS20171182.
2. Li F, et al. (2017). Apigenin Alleviates Endotoxin-Induced Myocardial Toxicity by Modulating Inflammation, Oxidative Stress, and Autophagy. *Oxid Med Cell Longev*. 2017:2302896. doi: 10.1155/2017/2302896.
3. Strickland, F.M. et al. (2017). Oxidative T cell modifications in lupus and Sjogren's syndrome. *Lupus* **2**(1):121.
4. Casós, K., et al (2017). Response of the human myocardium to ischemic injury and preconditioning: The role of cardiac and comorbid conditions, medical treatment, and basal redox status. *PLoS One*. **12**(4):e0174588. doi: 10.1371/journal.pone.0174588. eCollection 2017.
5. Lambrechts, K. et al. (2017). Venous gas emboli are involved in post-dive macro, but not microvascular dysfunction. *Eur J Appl Physiol*. **117**(2):335-344. doi: 10.1007/s00421-017-3537-9.
6. Shafique, E. et al. (2017). Mitochondrial redox plays a critical role in the paradoxical effects of NAPDH oxidase-derived ROS on coronary endothelium. *Cardiovasc Res*. **113**(2):234-246. doi: 10.1093/cvr/cvw249.
7. Wu P, et al. (2017). Mdivi-1 Alleviates Early Brain Injury After Experimental Subarachnoid Hemorrhage in Rats, Possibly via Inhibition of Drp1-Activated Mitochondrial Fission and Oxidative Stress. *Neurochem Res*. doi: 10.1007/s11064-017-2201-4.
8. Mikolka, P. et al. (2016). Lung inflammatory and oxidative alterations after exogenous surfactant therapy fortified with budesonide in rabbit model of meconium aspiration syndrome. *Physiol. Res*. **65** (Suppl. 5): S653-S662.
9. Ciarlone, G.E. and Dean, J.B. (2016). Acute hypercapnic hyperoxia stimulates reactive species production in the caudal solitary complex of rat brain slices but does not induce oxidative stress. *Am. J. Physiol. Cell Physiol*. **311**:C1027-C1039.
10. Mukhopadhyay, P. et al. (2016). PARP inhibition protects against alcoholic and nonalcoholic steatohepatitis. *J. Hepatol*. doi:10.1016/j.jhep.2016.10.023.
11. Bavunoglu, I. et al. (2016). Oxidative stress parameters and inflammatory and immune mediators as markers of the severity of sepsis. *J. Inf. Dev. Ctries*. **10**:1045-1052.
12. Liu, H. et al. (2016). Pterostilbene attenuates early brain injury following subarachnoid hemorrhage via inhibition of the NLRP3 inflammasome and Nox2-related oxidative stress. *Mol Neurobiol*. doi:10.1007/s12035-016-0108-8.
13. Saenen, N. D. et al. (2016). Lower placental leptin promoter methylation in association with fine particulate matter air pollution during pregnancy and placental nitrosative stress at birth in the ENVIRONAGE cohort. *Environ Health Perspect*. doi:10.1289/EHP38.

14. Martinez, P. F. et al. (2016). Modulation of MAPK and NF- κ B signaling pathways by antioxidant therapy in skeletal muscle of heart failure rats. *Cell Physiol Biochem*. doi:10.1159/000445631.
15. Dhall, S. et al. (2016). Tobacco toxins deposited on surfaces (third hand smoke) impair wound healing. *Clin Sci*. doi:10.1042/CS20160236.
16. Mikolka, P. et al. (2016). Antiinflammatory effect of N-Acetylcysteine combined with exogenous surfactant in meconium-induced lung injury. doi:10.1007/5584_2016_15.
17. Chan, T. K. et al. (2016). House dust mite-induced asthma causes oxidative damage and DNA double-strand breaks in the lungs. *J Allergy Clin Immunol*. doi:10.1016/j.jaci.2016.02.017.
18. Li, F. et al. (2016). Role of TFEB mediated autophagy, oxidative stress, inflammation, and cell death in endotoxin induced myocardial toxicity of young and aged mice. *Oxid Med Cell Longev*. doi:10.1155/2016/5380319.
19. Paul, S. et al. (2016). Size restricted Silymarin suspension evokes integrated adaptive response against acute hypoxia exposure in rat lung. *Free Radic Biol Med*. doi:10.1016/j.freeradbiomed.2016.04.020.
20. Adhami, N. et al. (2016). A health threat to bystanders living in the homes of smokers: How smoke toxins deposited on surfaces can cause insulin resistance. *PLoS One*. **11**:e0149510.
21. Unal, Betul, et al. (2016). Inhibition of neutral sphingomyelinase decreases elevated levels of nitrative and oxidative stress markers in liver ischemia-reperfusion injury. *Redox Rep*. doi:10.1080/13510002.2016.1162431.
22. Sumayao, R. et al. (2016). Lysosomal cystine accumulation promotes mitochondrial depolarization and induction of redox-sensitive genes in human kidney proximal tubular cells. *J Physiol*. doi:10.1113/JP271858.
23. Zhang, Z. Y. et al. (2015). Enhanced therapeutic potential of nano-curcumin against subarachnoid hemorrhage-induced blood-brain barrier disruption through inhibition of inflammatory response and oxidative stress. *Mol Neurobiol*. doi:10.1007/s12035-015-9635-y.

Please see the complete list of product citations: <http://www.cellbiolabs.com/nitrotyrosine-elisa-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

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