

---

Product Manual

# OxiSelect™ Total Glutathione (GSSG/GSH) Assay Kit

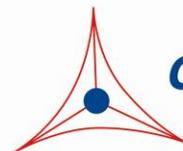
Catalog Number

STA-312

100 assays

**FOR RESEARCH USE ONLY**  
**Not for use in diagnostic procedures**

---



**CELL BIOLABS, INC.**  
*Creating Solutions for Life Science Research*

## **Introduction**

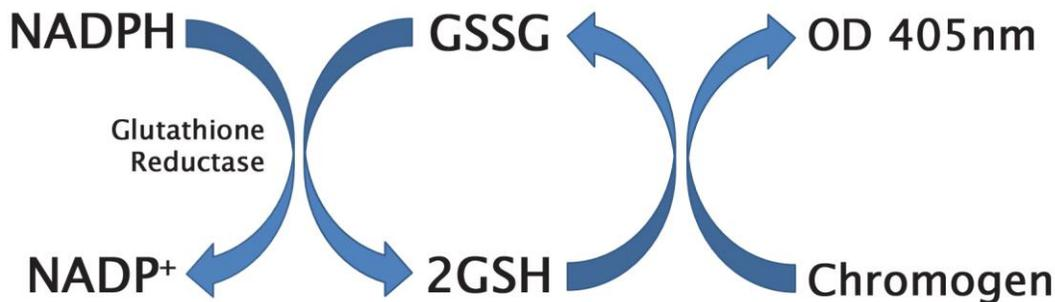
Glutathione is a key intracellular tripeptide thiol composed of glutamic acid, cysteine, and glycine. Glutathione helps protect cells from free radical damage by acting as an antioxidant. Within cells, glutathione exists in reduced (GSH) and oxidized (GSSG) states. In healthy cells and tissue, more than 90% of the total glutathione pool is in the reduced form (GSH) while less than 10% exists in the disulfide form (GSSG). The high GSH concentration is because the enzyme that transitions it from its oxidized state (GSSG), glutathione reductase, is constitutively active and inducible upon oxidative stress. An increased GSSG-to-GSH ratio is considered indicative of oxidative stress. Reduced glutathione's thiol group provides reducing equivalents to other unstable ROS, which in turn becomes unstable itself. This unstable GSH readily reacts with another unstable GSH to form a stable GSSG molecule. This reaction is prevalent since glutathione is present in high concentrations. GSSG is subsequently converted to GSH again by the enzyme glutathione reductase.

In addition to its role in oxidative stress, glutathione also helps maintain exogenous antioxidants such as vitamins C and E. Glutathione is involved with the breakdown of peroxides. It has a role in regulating the nitric acid cycle. Glutathione can directly bind many inorganic and organic xenobiotic (foreign chemicals) and carcinogenic compounds, such as the heavy metals mercury and arsenic. It is important to the proper function and maximum effect of the immune system. In addition, glutathione is fundamentally involved with many metabolic and biochemical mechanisms such as protein and prostaglandin synthesis, DNA synthesis and repair, maintenance of disulfide bonds in proteins, enzyme activation and amino acid transport across cell membranes.

Cell Biolabs' OxiSelect™ Total Glutathione Assay Kit is a quantitative assay for measuring the total concentration of glutathione, which encompasses both reduced and oxidized glutathione (GSH and GSSG) from plasma, blood, saliva, urine, tissue extracts, and plant or mammalian cell lysates. The kit employs a simple enzymatic recycling reaction for total glutathione quantification. The kit has a detection sensitivity limit of 8 nM. Each kit provides sufficient reagents to perform up to 100 assays, including standard curve and unknown samples.

## **Assay Principle**

The OxiSelect™ Total Glutathione Assay Kit is a quantitative assay for measuring the total glutathione content within a sample (GSH/GSSG). Glutathione Reductase reduces oxidized glutathione (GSSG) to reduced glutathione (GSH) in the presence of NADPH. Subsequently, the chromogen reacts with the thiol group of GSH to produce a colored compound that absorbs at 405 nm (Figure 1). The total glutathione content in unknown samples is determined by comparison with the predetermined glutathione standard curve. The rate of chromophore production is proportional to the concentration of glutathione within the sample. The rate can be determined from the absorbance change over time. Metaphosphoric acid is provided to remove interfering proteins or enzymes from samples.



**Figure 1. Assay Principle**

### **Related Products**

1. MET-5070: Glycine Assay Kit (Fluorometric)
2. MET-5054: L-Amino Acid Assay Kit (Colorimetric)
3. MET-5136: D-Amino Acid Assay Kit (Colorimetric)
4. MET-5080: Glutamate Assay Kit (Colorimetric)
5. STA-812: OxiSelect™ Glutathione Reductase Assay Kit (Colorimetric)

### **Kit Components**

#### **Box 1 (shipped at room temperature)**

1. Glutathione Reductase (50X) (Part No. 231201): One 50  $\mu$ L amber tube.
2. Chromogen (15X) (Part No. 231202): One 1 mL amber tube.
3. Assay Buffer (5X) (Part No. 231204): One 50 mL bottle.
4. Metaphosphoric Acid (MPA) (Part No. 231205): One 2 g bottle of crystals.
5. Glutathione Disulfide (GSSG) (Part No. 231206): One 50  $\mu$ L amber tube of a 1 mM solution.

#### **Box 2 (shipped on blue ice packs)**

1. NADPH (50X) (Part No. 231203): One 50  $\mu$ L amber tube.

### **Materials Not Supplied**

1. 96-well microtiter plate
2. Distilled or deionized water
3. 1X PBS

### **Storage**

Upon receipt, store the NADPH at -80°C. Prepare single use aliquots and avoid multiple freeze/thaw cycles. Store the remaining kit components at 4°C.

## **Preparation of Reagents**

- 1X Assay Buffer: Prepare 1X Assay Buffer by adding 200 mL of deionized water to 50 mL of the 5X Assay Buffer. Mix thoroughly until homogeneous. Use this buffer for preparing kit reagents. Store at 4°C when not in use.
- 1X Glutathione Reductase: Prepare the 1X Glutathione Reductase by diluting the stock solution 1:50 with 1X Assay Buffer. Vortex the stock tube thoroughly prior to preparing. Prepare only enough for immediate applications (eg. Add 25 µL of Glutathione Reductase stock to 1.225 mL Assay Buffer).
- 1X Chromogen: Prepare the 1X Chromogen just before use. Prepare only enough for immediate applications. Dilute the Chromogen stock 1:15 with 1X Assay Buffer (eg. Add 167 µL of Chromogen stock to 2.333 mL of 1X Assay Buffer. Vortex thoroughly.
- 1X NADPH: Prepare 1X NADPH by diluting the stock solution 1:50 with 1X Assay Buffer. Vortex the stock tube thoroughly prior to preparing. Prepare only enough for immediate applications (eg. Add 25 µL of NADPH stock to 1.225 mL Assay Buffer).
- Metaphosphoric Acid (MPA): Prepare a 5% (w/v) Metaphosphoric Acid (MPA) solution in deionized water. Prepare just before use. Prepare only enough for immediate applications (eg. Add 0.5 g of Metaphosphoric Acid crystals to 10 mL of deionized water). Vortex thoroughly.

*Note: MPA is corrosive and may cause burns. Use caution when handling acidic reagents.*

## **Preparation of Samples**

These preparation protocols are intended as a guide for preparing known samples. The user may need to adjust how they treat their sample accordingly. All samples should be assayed immediately or store at -80°C for up to 1-2 months. A trial assay with a representative test sample should be assayed to determine the samples compatibility with the dynamic range of the standard. It is recommended that samples be processed as soon as possible because GSH is rapidly metabolized and will continue to form various disulfides. The assay can be used on cell culture supernatants and lysates, blood, plasma, urine, saliva, tissue homogenates as well as other biological fluids. All samples should be treated with 5% MPA to remove interfering proteins and enzymes. The MPA treated deproteinated samples improve the stability of GSH. High levels of interfering substances may cause variations in results. Run proper controls as necessary. Always run a standard curve with samples.

### *Notes:*

- *Thiol compounds, such as cysteine, dithiothreitol (DTT), or β-mercaptoethanol can interfere with the assay by competing with GSH for binding to the Chromogen. In addition, N-ethylmaleimide or other thiol alkylating reagents should also be avoided because they will interfere with Glutathione Reductase and GSH.*
- *5% MPA will interfere with the assay. Upon preparing samples, each sample must be diluted 1:10 with 1X Assay Buffer to bring the MPA concentration to 0.5% final. Some samples may need to be diluted more. Make serial dilutions of samples as necessary to obtain a quantifiable change in absorbance readings over time.*
- *A kinetic assay is recommended because it is more precise than an end point assay.*
- **Saliva, Plasma or Urine:** GSSG in normal resting saliva, plasma and urine is at or below the detection limit for most glutathione assays. Collect sample in a microfuge tube and immediately add 4 volumes of ice-cold 5% MPA. Mix thoroughly and store on ice for 15 minutes. Centrifuge at 12,000 rpm for 10 minutes at 4°C to remove insoluble particles. Collect the supernatant. Store on ice if assaying immediately or freeze at -80°C for future use.

- **Cell Lysate:** Detach adherent cells by trypsinization. Count cells and centrifuge at 500 rpm for 5 minutes at 4°C. Wash cells with cold 1X PBS. Centrifuge suspension cells at 500 rpm for 5 minutes at 4°C. Remove supernatant and wash cells with cold 1X PBS. Repeat centrifugation and remove solution. Immediately resuspend the pellet with 200-500  $\mu$ L ice-cold 5% MPA for a cell concentration of  $1-5 \times 10^6$  cells. Mix thoroughly. Homogenize or sonicate cell suspension and store on ice until use. Transfer the suspension to a microfuge tube and centrifuge at 12,000 rpm for 5 minutes at 4°C. Collect the supernatant. Store on ice if used immediately or freeze at -80°C for future use.
- **Tissue Lysate:** The GSH concentration in most tissue is in the 1-10 mM range. It is recommended that a 10% w/v homogenate be created. Blood can contaminate a tissue sample due to high GSH concentrations. Therefore, perfusion of the tissue with a PBS/heparin is recommended. Remove tissue and wash the tissue thoroughly with cold isotonic saline solution of 1X PBS with 0.16 mg/mL heparin to prevent coagulation. Blot the tissue dry and weigh. Add ice-cold 5% MPA (~1 mL/100 mg tissue) and homogenize using a glass pestle. Centrifuge the homogenate at 12,000 rpm for 15 minutes at 4°C. Collect the supernatant. Store on ice if used immediately or freeze at -80°C for future use.
- **Erythrocyte Lysate:** Add blood sample to a blood collection tube with an anticoagulant such as heparin or sodium citrate. Centrifuge at 3,000 rpm for 15 minutes at 4°C. Remove and discard any plasma supernatant. Remove the white buffy coat (leukocytes) on the surface of the erythrocytes. Add four volumes of ice-cold 5% MPA to the pellet and resuspend. Mix thoroughly. Store on ice for 10 minutes. Centrifuge the suspension at 12,000 rpm for 10 minutes at 4°C. Collect the supernatant. Store on ice if used immediately or freeze at -80°C for future use.
- **Whole Blood Lysate:** Collect blood in conical tubes containing an anticoagulant such as sodium citrate or heparin. Add four volumes of ice-cold 5% MPA and mix thoroughly. Store on ice for 10 minutes. Centrifuge the suspension at 12,000 rpm for 10 minutes at 4°C. Collect the supernatant. Store on ice if used immediately or freeze at -80°C for future use.

### **Preparation of Standard Curve**

1. Prepare a solution of 0.5% MPA in 1X Assay Buffer and mix. Use this solution for preparing the glutathione standards.
2. To prepare the glutathione standards, first perform a 1:1000 dilution of the stock Glutathione Disulfide (GSSG) in 0.5% MPA/Assay Buffer. Use only enough for immediate applications (eg. Add 2  $\mu$ L of Glutathione Disulfide to 1998  $\mu$ L 0.5% MPA /Assay Buffer). This solution has a concentration of 1  $\mu$ M.
3. Use microfuge tubes to prepare a series of standards according to Table 1 below. Prepare standards fresh for each assay performed. Do not store or reuse standard preparations.

<b>Standard Tubes</b>	<b>1 <math>\mu</math>M GSSG Standard (<math>\mu</math>L)</b>	<b>0.5% MPA in Assay Buffer (<math>\mu</math>L)</b>	<b>GSSG (<math>\mu</math>M)</b>
1	500	500	0.5
2	500 of Tube #1	500	0.25
3	500 of Tube #2	500	0.125
4	500 of Tube #3	500	0.0625
5	500 of Tube #4	500	0.03125
6	500 of Tube #5	500	0.0156
7	500 of Tube #6	500	0.0078
8	0	500	0

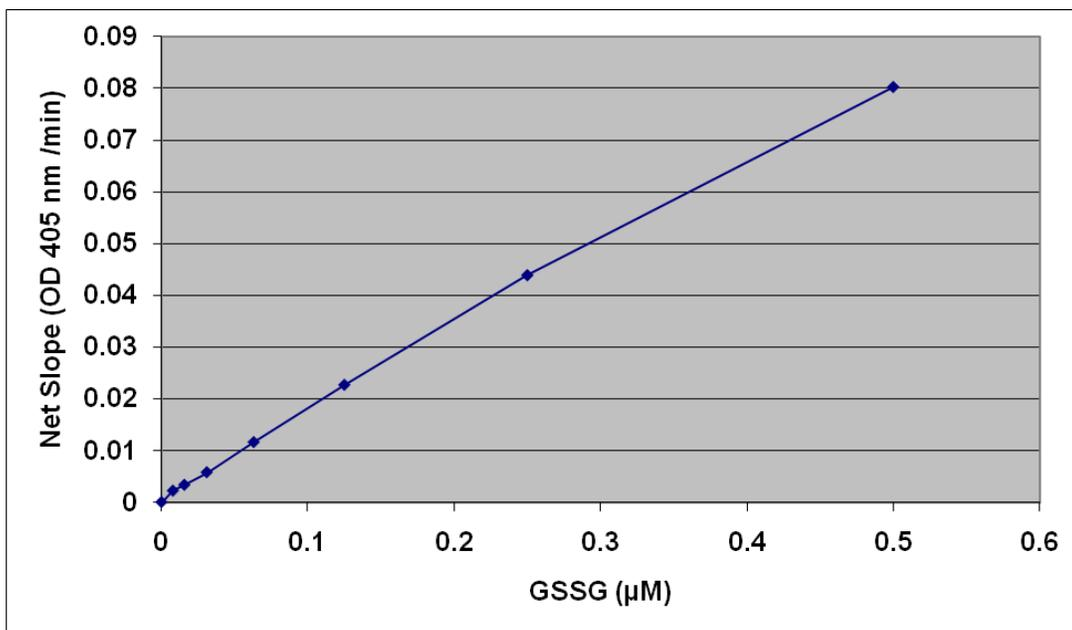
**Table 1. Preparation of Glutathione Disulfide Standards**

## **Total Glutathione (GSSG/GSH) Assay Protocol**

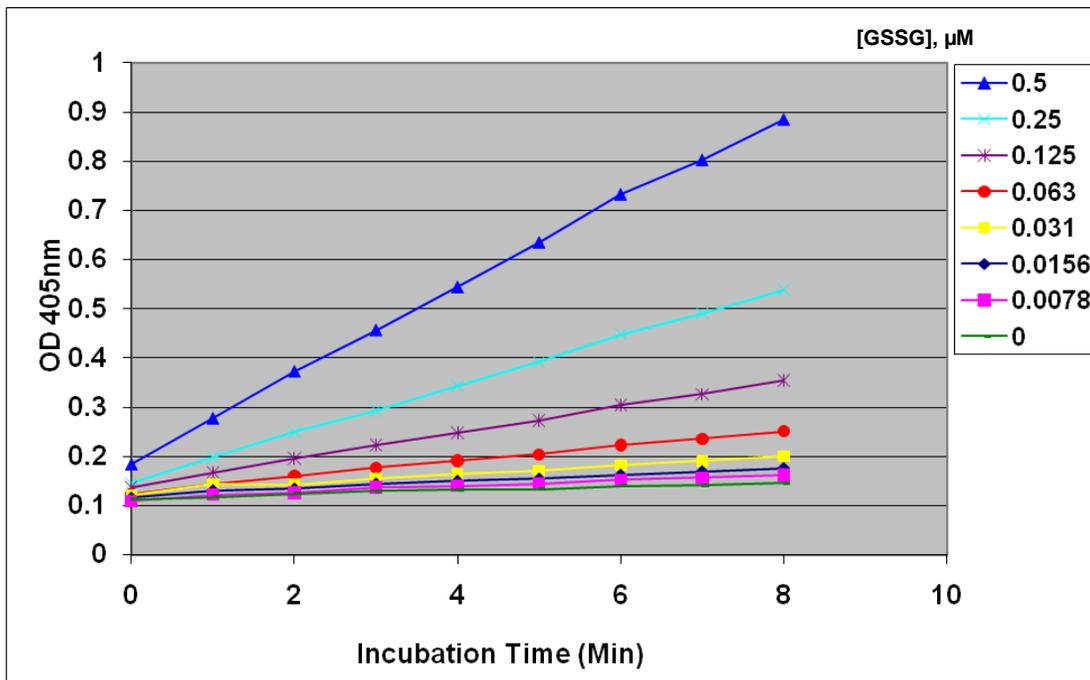
1. Prepare and mix all reagents thoroughly before use. Prepare the glutathione standards simultaneously with the samples so they may be assayed together. Each sample, including unknown and standard, should be assayed in duplicate or triplicate.
2. In a 96-well plate, add 25  $\mu\text{L}$  of the 1X Glutathione Reductase solution to each well to be tested.
3. Add 25  $\mu\text{L}$  of the 1X NADPH solution to each well to be tested.
4. Add 100  $\mu\text{L}$  of the prepared glutathione standards or samples to each well to be tested. Mix thoroughly.
5. Ensure that the plate reader is prepared for a kinetic assay and is set to read at 405 nm.
6. Add 50  $\mu\text{L}$  of the 1X Chromogen and mix briefly. Immediately begin recording the absorbance at 405 nm at 1-minute intervals for 10 minutes. If using all the wells within the plate at one time, then it may be necessary to record the absorbance at 2-minute intervals.
7. Calculate the concentration of standards and samples. See Calculation of Results below.

## **Example of Results**

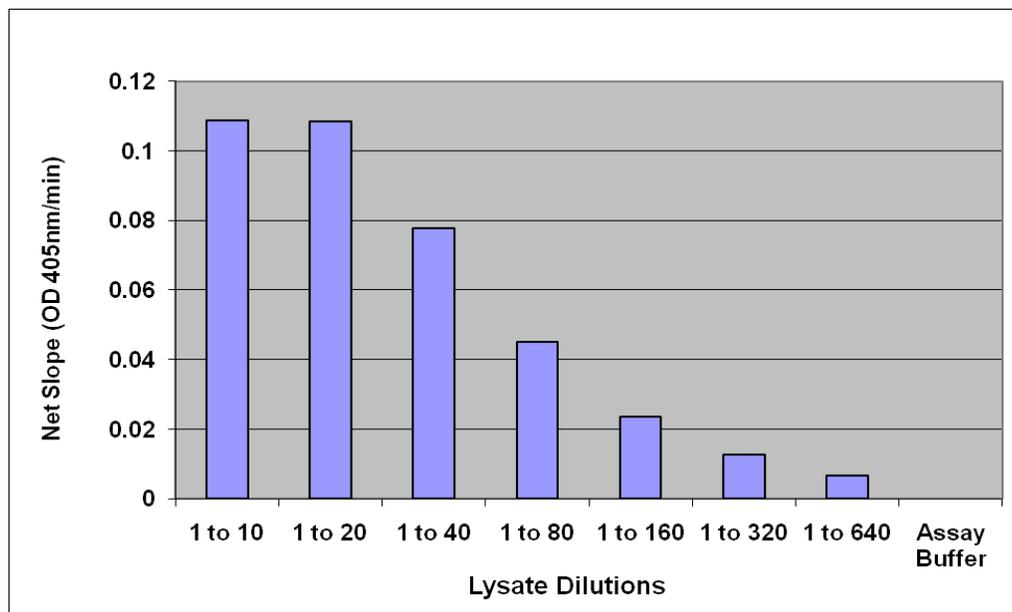
The following figures demonstrate typical Total Glutathione Assay results. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 3: GSSG Standard Curve (Net slope versus GSSG concentration).**



**Figure 2. GSSG Standard Curve** (OD 405nm versus incubation time as a function of GSSG concentration).



**Figure 4. Total Glutathione Content in NIH 3T3 Cell Lysate.** Samples were prepared according to the instructions under Sample Preparations above.

### Calculation of Results

1. First, determine the average of the replicate absorbance readings for each GSSG standard, sample, and negative control for every time point taken.
2. Graph the average of each standard, sample, and background absorbance at 405nm against incubation time. Determine the slope for each value from the linear portion of each curve. See Figure 2 under Example of Results below.
3. Next, subtract the background slope from the slope of the standards and samples.

4. Plot the net slopes of the GSSG standards against the micromolar concentration of GSSG. See Figure 3 under Example of Results below.
5. Compare the net slopes of the samples with the standard curve from Figure 3 and determine the micromolar concentration of GSSG for each sample. This is equivalent to the total glutathione content.

## **References**

1. Anderson, M. Glutathione in Free radicals, A Practical Approach. Oxford University Press, New York; 1996.
2. Halliwell, B, Gutteridge, J.M.C. Free Radicals in Biology and Medicine. Oxford University Press, New York; 1999.
3. Julius, M., et al. *J. Clin. Epidemiol.* (1994) 47: 1021-1026.
4. Mytilineou, C., et al. *Parkinsonism Relat. Disord.* (2002) 8: 385-387.

## **Recent Product Citations**

1. Al-Masri, A. A. et al. (2023). Anti-inflammatory effect of anthocyanin-rich extract from banana bract on lipopolysaccharide-stimulated RAW 264.7 macrophages. *J Funct Foods*. doi: 10.1016/j.jff.2023.105628.
2. Klyosova, E. et al. (2022). A Polymorphism in the Gene Encoding Heat Shock Factor 1 (HSF1) Increases the Risk of Type 2 Diabetes: A Pilot Study Supports a Role for Impaired Protein Folding in Disease Pathogenesis. *Life (Basel)*. **12**(11):1936. doi: 10.3390/life12111936.
3. Torres-Vargas, J.A. et al. (2022). Characterization of the activity and the mechanism of action of a new toluquinol derivative with improved potential as an antiangiogenic drug. *Biomed Pharmacother*. doi: 10.1016/j.biopha.2022.113759.
4. Althobaiti, N.A. et al. (2022). Ethnomedicinal Plants with Protective Effects against Beta-Amyloid Peptide (A $\beta$ )1-42 Indicate Therapeutic Potential in a New In Vivo Model of Alzheimer's Disease. *Antioxidants*. **11**(10):1865. doi: 10.3390/antiox11101865.
5. Rajab, B.S. et al. (2022). Antioxidative and Anti-Inflammatory Protective Effects of  $\beta$ -Caryophyllene against Amikacin-Induced Nephrotoxicity in Rat by Regulating the Nrf2/AMPK/AKT and NF- $\kappa$ B/TGF- $\beta$ /KIM-1 Molecular Pathways. *Oxid Med Cell Longev*. doi: 10.1155/2022/4212331.
6. Tsunenaga, M. et al. (2022). Modulating effects of oral administration of Lycii Fructus extracts on UVB-induced skin erythema: A Randomized, placebo-controlled study. *Biomed Rep*. **17**(1):62. doi: 10.3892/br.2022.1545.
7. Polonikov, A. et al. (2022). The Impact of Genetic Polymorphisms in Glutamate-Cysteine Ligase, a Key Enzyme of Glutathione Biosynthesis, on Ischemic Stroke Risk and Brain Infarct Size. *Life (Basel)*. **12**(4):602. doi: 10.3390/life12040602.
8. Giri, T. et al. (2022). Labor induction with oxytocin in pregnant rats is not associated with oxidative stress in the fetal brain. *Sci Rep*. **12**(1):3143. doi: 10.1038/s41598-022-07236-x.
9. Refaat, B. & El-Boshy, M. (2022). Protective antioxidative and anti-inflammatory actions of  $\beta$ -caryophyllene against sulfasalazine-induced nephrotoxicity in rat. *Exp Biol Med (Maywood)*. doi: 10.1177/15353702211073804.
10. Sommerfeld-Klatta, K. et al. (2022). Oxidative stress and biochemical indicators in blood of patients addicted to alcohol treated for acute ethylene glycol poisoning. *Hum Exp Toxicol*. **41**:9603271211061502. doi: 10.1177/09603271211061502.
11. Olszewska-Czyz, I. et al. (2022). The Influence of Hyaluronic Acid Adjunctive Therapy of Periodontitis on Salivary Markers of Oxidative Stress: Randomized, Controlled Clinical Trial. *Antioxidants (Basel)*. **11**(1):135. doi: 10.3390/antiox11010135.

12. Zhang, Q. et al. (2021). Micelle silymarin supplementation to sows' diet from day 109 of gestation to entire lactation period enhances reproductive performance, and affects serum hormones and metabolites. *J Anim Sci*. doi: 10.1093/jas/skab354.
13. Kalinovic, S. et al. (2021). Comparison of three methods for in vivo quantification of glutathione in tissues of hypertensive rats. *Free Radic Res*. doi: 10.1080/10715762.2021.2016735.
14. Majewski, M. et al. (2021). The Role of 20-HETE, COX, Thromboxane Receptors, and Blood Plasma Antioxidant Status in Vascular Relaxation of Copper-Nanoparticle-Fed WKY Rats. *Nutrients*. **13**(11):3793. doi: 10.3390/nu13113793.
15. Park, H.S. et al. (2021). Gastroprotective Effect of Myricetin on Ethanol-Induced Acute Gastric Injury in Rats. *Evid Based Complement Alternat Med*. doi: 10.1155/2021/9968112.
16. Althobaiti, N.A. et al. (2021). Assessment and Validation of *Globodera pallida* as a Novel In Vivo Model for Studying Alzheimer's Disease. *Cells*. **10**(9):2481. doi: 10.3390/cells10092481.
17. 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.
18. Azarova, I. et al. (2021). The Link between Type 2 Diabetes Mellitus and the Polymorphisms of Glutathione-Metabolizing Genes Suggests a New Hypothesis Explaining Disease Initiation and Progression. *Life (Basel)*. **11**(9):886. doi: 10.3390/life11090886.
19. Kozłowski, K. et al. (2021). Growth performance, immune status and intestinal fermentative processes of young turkeys fed diet with additive of full fat meals from *Tenebrio molitor* and *Hermetia illucens*. *Anim Feed Sci Technol*. doi: 10.1016/j.anifeeds.2021.114994.
20. Almarhoun, M. et al. (2021). Overexpression of STARD3 attenuates oxidized LDL-induced oxidative stress and inflammation in retinal pigment epithelial cells. *Biochim Biophys Acta Mol Cell Biol Lipids*. **1866**(7):158927. doi: 10.1016/j.bbalip.2021.158927.
21. Gupta, S. et al. (2021). Glutathione is a potential therapeutic target for acrolein toxicity in the cornea. *Toxicol Lett*. doi: 10.1016/j.toxlet.2021.01.005.
22. Nour, M.A. et al. (2021). Productive performance, fertility and hatchability, blood indices and gut microbial load in laying quails as affected by two types of probiotic bacteria. *Saudi J Biol Sci*. doi: 10.1016/j.sjbs.2021.07.030.
23. Kikuchi, M. et al. (2021). Effects of glucoraphanin-rich broccoli sprout extracts on sleep quality in healthy adults: An exploratory study. *J Funct Foods*. doi: 10.1016/j.jff.2021.104574.
24. Jiang, J. et al. (2021). Impact of intrauterine fetal resuscitation with oxygen on oxidative stress in the developing rat brain. *Sci Rep*. **11**(1):9798. doi: 10.1038/s41598-021-89299-w.
25. 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.
26. de Los Santos-Jiménez, J. et al. (2021). Glutaminase isoforms expression switches microRNA levels and oxidative status in glioblastoma cells. *J Biomed Sci*. **28**(1):14. doi: 10.1186/s12929-021-00712-y.
27. Rafeeq, M. et al. (2021). Protective effect of 6-paradol in acetic acid-induced ulcerative colitis in rats. *BMC Complement Med Ther*. **21**(1):28. doi: 10.1186/s12906-021-03203-7.
28. Zapata-Londono, M.B. et al. (2020). Effect of mango (*Mangifera indica*) cv. azÚcar juice consumption on plasma antioxidant capacity and oxidative stress biomarkers. *Vitae*. **27**(1):1-10. doi: 10.17533/udea.vitae.v27n1a03.
29. Sanz-González, S.M. et al. (2020). Clinical and Molecular-Genetic Insights into the Role of Oxidative Stress in Diabetic Retinopathy: Antioxidant Strategies and Future Avenues. *Antioxidants (Basel)*. **9**(11): E1101. doi: 10.3390/antiox9111101.

30. Bakalova, R. et al. (2020). Selective Targeting of Cancerous Mitochondria and Suppression of Tumor Growth Using Redox-Active Treatment Adjuvant. *Oxid Med Cell Longev*. doi: 10.1155/2020/6212935.

### **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.  
5628 Copley Drive  
San Diego, CA 92111  
Worldwide: +1 858 271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2010-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.