
Product Manual

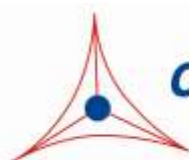
OxiSelect™ Total Antioxidant Capacity (TAC) Assay Kit

Catalog Number

STA-360

200 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

Oxidative stress is a physiological condition where there is an imbalance between concentrations of reactive oxygen species (ROS) and antioxidants. However, excessive ROS accumulation will lead to cellular injury, such as damage to DNA, proteins, and lipid membranes. The cellular damage caused by ROS has been implicated in the development of many disease states, such as cancer, diabetes, cardiovascular disease, atherosclerosis, and neurodegenerative diseases. Under normal physiological conditions, cellular ROS generation is counterbalanced by the action of cellular antioxidant enzymes, macro or micro molecules, as well as other redox molecules. Antioxidants also include both hydrophilic and lipophilic molecules for metabolizing ROS. These may be localized transiently within different tissues or cells. Because of their potential harmful effects, excessive ROS must be promptly eliminated from the cells by this variety of antioxidant defense mechanisms.

Although the products of ROS-induced oxidative stress are extensively used to monitor their biological effects, it is also important to evaluate the antioxidant capacity of biological fluids, cells, and extracts. Antioxidants commonly neutralize radicals via a hydrogen atom transfer (HAT) or single electron transfer (SET) mechanism. SET assays quantify the capability of an antioxidant to transfer one electron to reduce any compound, such as free radicals, carbonyls, and metals.

Cell Biolabs' OxiSelect™ Total Antioxidant Capacity (TAC) Assay Kit measures the total antioxidant capacity of biomolecules from a variety of samples via a SET mechanism. The TAC Assay is based on the reduction of copper (II) to copper (I) by antioxidants such as uric acid. Upon reduction, the copper (I) ion further reacts with a coupling chromogenic reagent that produces a color with a maximum absorbance at 490 nm. The net absorbance values of antioxidants are compared with a known uric acid standard curve. Absorbance values are proportional to the sample's total reductive capacity. Results are expressed as "µM Copper Reducing Equivalents" or "mM Uric Acid Equivalents". Copper is advantageous over iron-based antioxidant assays because all classes of antioxidants, including thiols, are detected with marginal radical interference. In addition, the kinetics of the copper-based reaction is also faster than the iron-based reaction, which makes the TAC Assay an ideal tool for estimating reductive or antioxidant capacity efficiently and accurately.

Cell Biolabs' OxiSelect™ TAC Assay Kit is a fast and reliable kit for the direct measurement of total antioxidant capacity from cell lysate, plasma, serum, urine, tissue homogenates, and food extracts. Each kit provides sufficient reagents to perform up to 200 assays, including blanks, antioxidant standards and unknown samples. Hydrophilic and lipophilic samples are compatible with the assay.

Assay Principle

Cell Biolabs' OxiSelect™ TAC Assay Kit measures the total antioxidant capacity within a sample. Samples are compared to a known concentration of uric acid standard within a 96-well microtiter plate format. Samples and standards are diluted with a reaction reagent and, upon the addition of copper, the reaction proceeds for a few minutes. The reaction is stopped and read with a standard 96-well spectrophotometric microplate reader at 490 nm (Figure 1). Antioxidant capacity is determined by comparison with the uric acid standards.

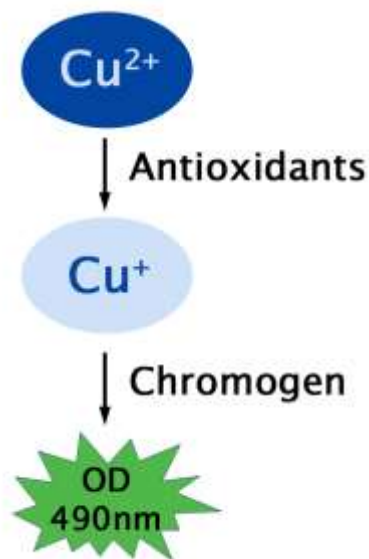


Figure 1. TAC Activity Assay Principle

Related Products

1. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit
2. STA-312: OxiSelect™ Total Glutathione (GSSG/GSH) Assay Kit
3. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
4. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
5. STA-340: OxiSelect™ Superoxide Dismutase Activity Assay
6. STA-341: OxiSelect™ Catalase Activity Assay
7. STA-342: OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence)
8. STA-344: OxiSelect™ Hydrogen Peroxide/Peroxidase Assay (Fluorometric)
9. STA-345: OxiSelect™ ORAC Activity Assay
10. STA-346: OxiSelect™ HORAC Activity Assay
11. STA-347: OxiSelect™ In Vitro ROS/RNS Assay Kit (Green Fluorescence)

Kit Components

1. Uric Acid Standard (Part No. 236001): One 100 mg tube of powder.
2. Reaction Buffer (100X) (Part No. 236002): One 400 μL amber tube.
3. Copper Ion Reagent (100X) (Part No. 236003): One 1.0 mL tube.
4. Stop Solution (10X) (Part No. 236004): One 1.5 mL tube.

Materials Not Supplied

1. Standard 96-well microtiter plates for use in microplate reader
2. 1N NaOH, 1X PBS and deionized water
3. Methanol or other organic solvent for lipid-based samples
4. Sonicator or homogenizer for sample preparations
5. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
6. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
7. Spectrophotometric microplate reader capable of reading 490 nm

Storage

Upon receipt store the Reaction Buffer (100X) at 4°C. Store all remaining kit components at room temperature.

Preparation of Reagents

Reagents may be prepared for either hydrophilic or lipophilic samples. Although many lipophilic samples are soluble upon dilution with 1X PBS, kit reagents and samples may be prepared in methanol to ensure complete solubility.

- 1X Reaction Buffer: Dilute the Reaction Buffer 1:100 with 1X PBS (hydrophilic) or with methanol (lipophilic). Mix to homogeneity. Store the 1X Reaction Buffer at 4°C up to three months.
- 1X Copper Ion Reagent: Dilute the Copper Ion Reagent 1:100 with deionized water (hydrophilic) or with methanol (lipophilic). Mix to homogeneity. Store the 1X Copper Ion Reagent at 4°C up to three months.
- 1X Stop Solution: Dilute the Stop Solution 1:10 with deionized water (hydrophilic) or with methanol (lipophilic). Mix to homogeneity. Store the 1X Stop Solution at 4°C up to three months.

Preparation of Samples

Samples should be stored at -80°C prior to performing the assay. Samples should be prepared at the discretion of the user. The following recommendations are only guidelines and may be altered to optimize or complement the user's experimental design.

Note: EDTA can interfere with the TAC assay and should not be present in any sample.

- Tissue Lysate: Sonicate or homogenize tissue sample on cold PBS and centrifuge at 10,000 x g for 10 minutes at 4°C. Aliquot the supernatant for storage at -80°C, protein determination and subsequent TAC assay.
- Cell Culture: Wash cells 3 times with cold PBS prior to lysis. Lyse cells with sonication or homogenation in cold PBS and centrifuge at 10,000 x g for 10 minutes at 4°C. Aliquot the supernatant for storage at -80°C, protein determination and subsequent TAC assay.
- Plasma: Collect blood with heparin and centrifuge at 4°C for 10 minutes. Remove the plasma and aliquot samples for testing.
- Urine: Test neat or diluted with PBS if appropriate.

- Lipophilic Fractions: Dissolve lipophilic samples in 100% methanol or acetone and then dilute in 50% methanol or acetone. Incubate the mixture for 1 hour at room temperature with mixing. Further dilute samples as necessary prior to testing.
- Food Samples: Results may vary depending on sample source and purification. Dilution and preparation of these samples is at the discretion of the user, but use the following guidelines:
 - Solid or High Protein Samples: Weigh solid sample and then homogenize after adding deionized water (1:2, w/v). Centrifuge the homogenate at 10,000 x g for 10 minutes at 4°C. Recover the supernatant which is the water-soluble fraction. The insoluble fraction (pulp) is further extracted by adding pure acetone (1:4, w(solid pulp)/v) and mixing at room temperature for 30-60 minutes. Centrifuge the extract/solid at 10,000 x g for 10 minutes at 4°C. Recover the acetone extract and dilute with PBS or water as necessary prior to running the assay. The TAC value is calculated by combining the results from the water-soluble fraction and the acetone extract from the pulp fraction.
 - Aqueous Samples: Centrifuge the sample at 10,000 x g for 10 minutes at 4°C to remove any particulates. Dilute the supernatant in PBS as necessary prior to running the assay. Certain liquids such as juice extracts may be tested without dilution.

Preparation of Uric Acid Standard Curve

1. Prepare fresh Uric Acid standards by weighing out the Uric Acid powder for a 10 mg/mL solution in 1N NaOH. This 10 mg/mL is equivalent to a concentration of 60 mM. Use the 60 mM Uric Acid solution to prepare a 2 mM solution of Uric Acid (eg. add 100 μ L of the 60 mM Uric Acid standard to 2.900 mL of deionized water).
2. Prepare a series of the remaining Uric Acid standards according to Table 1 below. Prepare in deionized water for hydrophilic testing or methanol for lipophilic testing.

Tubes	2 mM Uric Acid Antioxidant Standard (μL)	Deionized Water or Methanol (μL)	Resulting Uric Acid Concentration (mM)
1	500	500	1
2	500 of tube #1	500	0.5
3	500 of tube #2	500	0.25
4	500 of tube #3	500	0.125
5	500 of tube #4	500	0.0625
6	500 of tube #5	500	0.03125
7	500 of tube #6	500	0.0156
8	500 of tube #7	500	0.0078
9	500 of tube #8	500	0.0039
10	0	500	0.0

Table 1. Preparation of Uric Acid Standards.

Note: The 60 mM stock Uric Acid standard solution may be aliquotted and stored at -70°C for up to one week. Do not store diluted Uric Acid Standard solutions.

Assay Protocol

Each Uric Acid Standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

1. Add 20 μL of the diluted Uric Acid Standards or samples to the 96-well microtiter plate.
2. Add 180 μL of the 1X Reaction Buffer to each well using either a multichannel pipette or a plate reader liquid handling system. Mix thoroughly.
3. Obtain an initial absorbance by reading the plate at 490 nm.
4. To initiate the reaction, add 50 μL of the 1X Copper Ion Reagent into each well. Incubate 5 minutes on an orbital shaker.
5. Add 50 μL of 1X Stop Solution to each well to terminate the reaction.
6. Read the plate again at 490 nm.

Example of Results

The following figures demonstrate typical OxiSelect™ TAC Assay results (hydrophilic). One should use the data below for reference only. This data should not be used to interpret or calculate actual sample results.

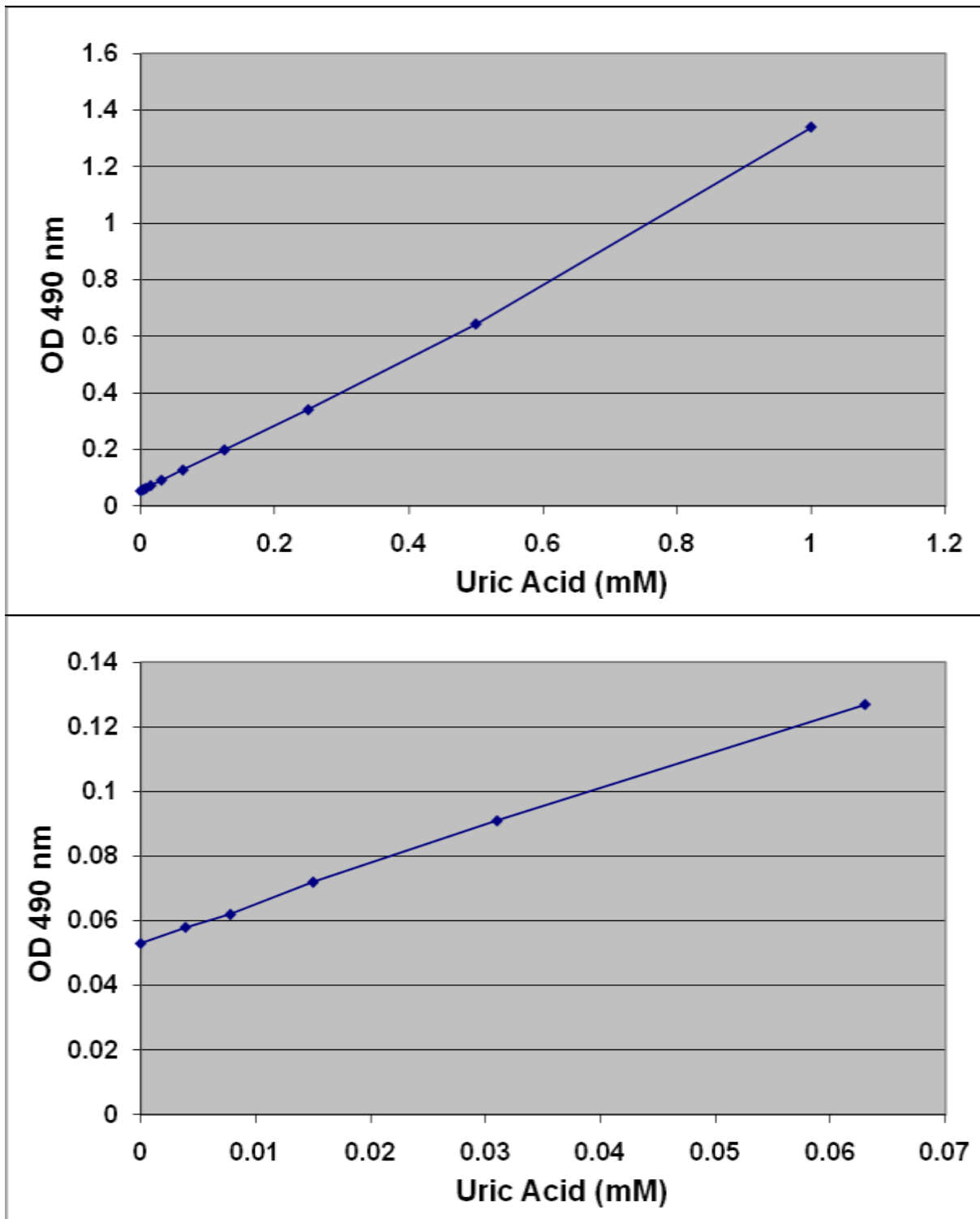


Figure 2: TAC Assay Standard Curve.

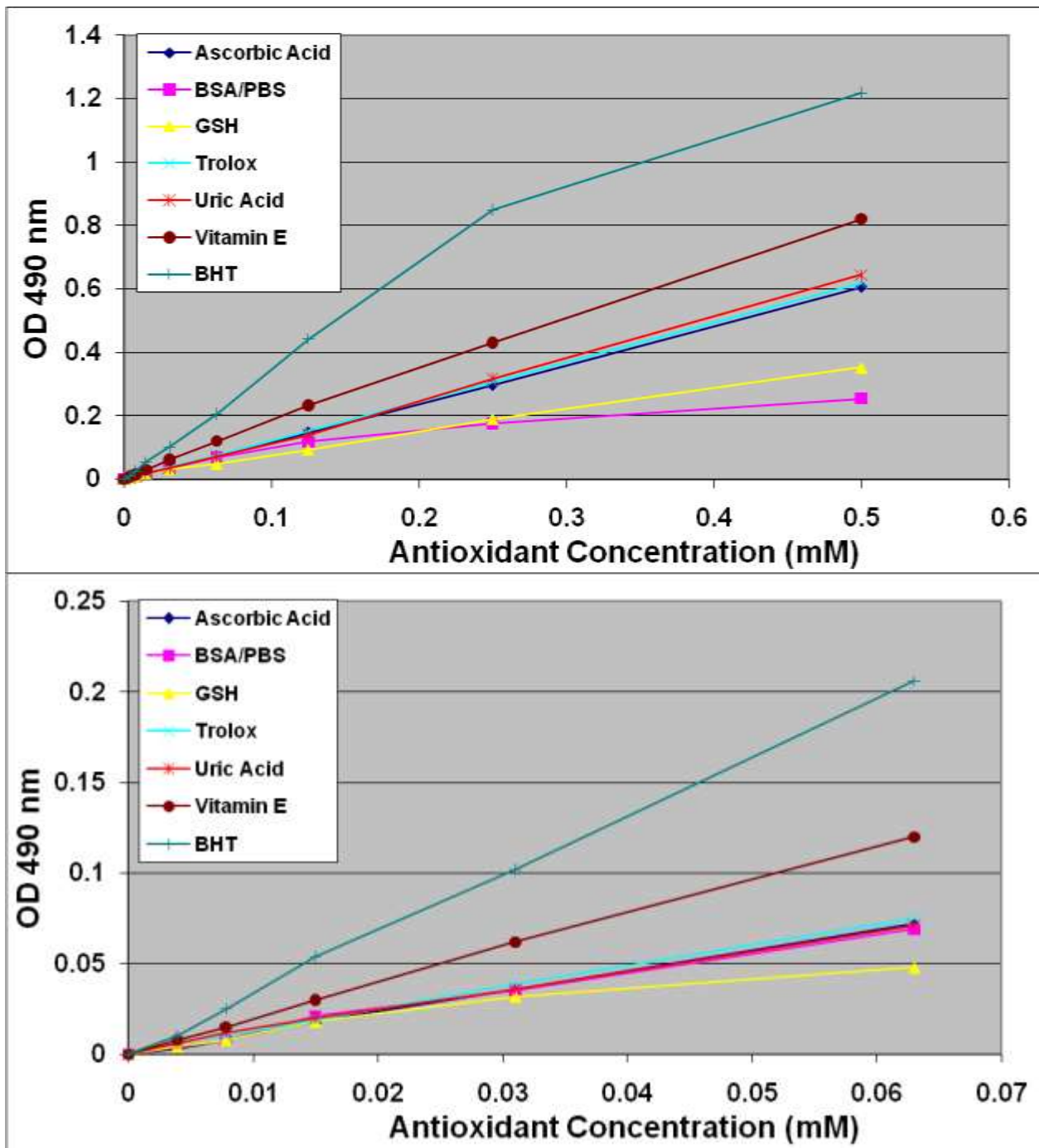


Figure 3: TAC Assay of various antioxidants.

Calculation of Results

1. Calculate the net absorbance by subtracting the initial absorbance readings for samples and standards (Step 3) from the final readings taken for each (Step 6).
2. Plot the net absorbance against the uric acid concentration for the uric acid standard curve.
3. Calculate the antioxidant capacity of unknown samples by comparing the net OD 490 nm values of samples to the uric acid standard curve.

To determine the “mM uric acid equivalents” (UAE) for samples, extrapolate the uric acid concentration from the sample’s analogous uric acid OD 490 nm value (For example, based on the sample standard curve above, an OD 490 nm value of 1.0 corresponds to 0.75 mM uric acid equivalents (UAE)).

To determine “ μ M Copper Reducing Equivalents” (CRE) for samples, multiply the uric acid equivalence (UAE) concentration by 2189 μ M Cu^{++} / mM uric acid. **1 mM of uric acid = 2189 μ M Copper Reducing Equivalents.** (For example, 0.75 mM UAE x 2189 = 1642 Copper Reducing Equivalents (CRE)). CRE sample values are proportional to the sample’s Total Antioxidant Capacity or Total Antioxidant Power.

References

1. Allard, J.P., et al. (1998) *Am. J. Clin. Nutr.* **67**: 143-147.
2. Cerutti, P. and Trump, B. (1991) *Proc. Cancer Cell* **3**: 1-7.
3. Frei, B., et al. (1992) “Molecular Biology of Free Radical Scavenging System” 23-45.
4. Trachootham, D., et al. (2008) *Antioxid. Redox Signal.* **10**: 1343-1374.
5. Van-Zoeren-Grobbe, et al. (1997) *Acta Paediatrica.* **86**: 1356-1362.

Recent Product Citations

1. Lingens, J.B. et al. (2021). Evaluation of Methionine Sources in Protein Reduced Diets for Turkeys in the Late Finishing Period Regarding Performance, Footpad Health and Liver Health. *Agriculture*. **11**(9):901. doi: 10.3390/agriculture11090901.
2. Ahmed-Farid, O.A. et al. (2021). Effects of Chronic Thermal Stress on Performance, Energy Metabolism, Antioxidant Activity, Brain Serotonin, and Blood Biochemical Indices of Broiler Chickens. *Animals (Basel)*. **11**(9):2554. doi: 10.3390/ani11092554.
3. Saber, T.M. et al. (2021). Early postmortem biochemical, histological, and immunohistochemical alterations in skeletal muscles of rats exposed to boldenone undecylenate: Forensic implication. *J Forensic Leg Med.* **83**:102248. doi: 10.1016/j.jflm.2021.102248.
4. Tembo, M.C. et al. (2021). The Predictability of Frailty Associated with Musculoskeletal Deficits: A Longitudinal Study. *Calcif Tissue Int.* doi: 10.1007/s00223-021-00865-w.
5. Delos Reyes, J.B. et al. (2021). Effects of dietary supplementation of vitamin C on productive performance, egg quality, tibia characteristics and antioxidant status of laying hens. *Livest Sci.* doi: 10.1016/j.livsci.2021.104502.
6. Tembo, M.C. et al. (2021). Association between serum interleukin-6 and frailty in older men: cross-sectional data. *Eur Geriatr Med.* doi: 10.1007/s41999-021-00490-8.
7. Abo El Gheit, R.E. et al. (2021). Retinoprotective effect of agmatine in streptozotocin-induced diabetic rat model: avenues for vascular and neuronal protection : Agmatine in diabetic retinopathy. *J Physiol Biochem.* doi: 10.1007/s13105-021-00799-9.

8. Serrya, M.S. et al. (2021). Evaluation of the therapeutic effects of mycophenolate mofetil targeting Nrf-2 and NLRP3 inflammasome in acetic acid induced ulcerative colitis in rats. *Life Sci.* doi: 10.1016/j.lfs.2021.119154.
9. Akgül, B. et al. (2021). Alleviation of prilocaine-induced epileptiform activity and cardiotoxicity by thymoquinone. *Daru.* doi: 10.1007/s40199-020-00385-2.
10. Yu, D.G. et al. (2021). Effects of stocking density and dietary vitamin C on performance, meat quality, intestinal permeability, and stress indicators in broiler chickens. *J Anim Sci Technol.* doi: 10.5187/jast.2021.e77.
11. Gross, J.J. et al. (2021). Metabolic and endocrine responses to short-term nutrient imbalances in the feed ration of mid-lactation dairy cows. *Animal.* **15**(7):100306. doi: 10.1016/j.animal.2021.100306.
12. Keller, M. et al. (2021). Antioxidant and Inflammatory Gene Expression Profiles of Bovine Peripheral Blood Mononuclear Cells in Response to *Arthrospira platensis* before and after LPS Challenge. *Antioxidants (Basel).* **10**(5):814. doi: 10.3390/antiox10050814.
13. El-Sadek, H.M. et al. (2021). Pentoxifylline treatment alleviates kidney ischemia/reperfusion injury: Novel involvement of galectin-3 and ASK-1/JNK & ERK1/2/NF- κ B/HMGB-1 trajectories. *J Pharmacol Sci.* doi: 10.1016/j.jphs.2021.03.011.
14. Bakeer, M.R. et al. (2021). Effect of dietary pumpkin (*Cucurbita moschata*) seed oil supplementation on reproductive performance and serum antioxidant capacity in male and nulliparous female V-Line rabbits. *Ital. J. Anim. Sci.* **20**(1):419-425. doi: 10.1080/1828051X.2021.1889406.
15. Lewandowski, L. et al. (2021). Concentration/activity of superoxide dismutase isozymes and the pro-/antioxidative status, in context of type 2 diabetes and selected single nucleotide polymorphisms (genes: INS, SOD1, SOD2, SOD3) – Preliminary findings. *Biomed Pharmacother.* doi: doi.org/10.1016/j.biopha.2021.111396.
16. 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.
17. Ra, K. et al. (2021). Comparison of Anti-Oxidative Effect of Human Adipose- and Amniotic Membrane-Derived Mesenchymal Stem Cell Conditioned Medium on Mouse Preimplantation Embryo Development. *Antioxidants (Basel).* **10**(2):268. doi: 10.3390/antiox10020268.
18. Satpathy, S. et al. (2021). A fraction of *Pueraria tuberosa* extract, rich in antioxidant compounds, alleviates ovariectomized-induced osteoporosis in rats and inhibits growth of breast and ovarian cancer cells. *PLoS One.* **16**(1):e0240068. doi: 10.1371/journal.pone.0240068.
19. 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.
20. Tsvetkova, A.S. et al. (2020). Melatonin Prevents Early but Not Delayed Ventricular Fibrillation in the Experimental Porcine Model of Acute Ischemia. *Int J Mol Sci.* **22**(1):E328. doi: 10.3390/ijms22010328.
21. Rondanelli, M. et al. (2020). Short- and Long-Term Effectiveness of Supplementation with Non-Animal Chondroitin Sulphate on Inflammation, Oxidative Stress and Functional Status in Obese Subjects with Moderate Knee Osteoarthritis before and after Physical Stress: A Randomized, Double-Blind, Placebo-Controlled Trial. *Antioxidants (Basel).* **9**(12):1241. doi: 10.3390/antiox9121241.
22. Domingo, J.C. et al. (2020). Are Circulating FGF21 and NT-proBNP promising novel biomarkers in Myalgic Encephalomyelitis/Chronic Fatigue Syndrome? *Antioxid Redox Signal.* doi: 10.1089/ars.2020.8230.

23. Manzocchi, E. et al. (2020). Effects of the substitution of soybean meal by spirulina in a hay-based diet for dairy cows on milk composition and sensory perception. *J Dairy Sci.* doi: 10.3168/jds.2020-18602.
24. Silva-Guillen, Y. et al. (2020). Growth performance, oxidative stress, and antioxidant capacity of newly weaned piglets fed dietary peroxidized lipids with vitamin E or phytochemical compounds in drinking water. *Applied Animal Science.* doi: 10.15232/aas.2019-01976.
25. Drobic, F. et al. (2020). Efficient Muscle Distribution Reflects the Positive Influence of Coenzyme Q10 Phytosome in Healthy Aging Athletes after Stressing Exercise. *J Food Sci Nutr Res.* **3**(4):262-275. doi: 10.26502/jfsnr.2642-11000054.
26. Moustafa, A. (2020). Effect of Light-Dark Cycle Misalignment on the Hypothalamic-Pituitary-Gonadal Axis, Testicular Oxidative Stress, and Expression of Clock Genes in Adult Male Rats. *Int J Endocrinol.* doi: 10.1155/2020/1426846.
27. Sen, U. et al. (2020). Differential sensitivities of triple-negative breast cancer stem cell towards various doses of vitamin C: An insight into the internal antioxidant systems. *J Cell Biochem.* doi: 10.1002/jcb.29863.
28. Guedj, F. et al. (2020). Apigenin as a Candidate Prenatal Treatment for Trisomy 21: Effects in Human Amniocytes and the Ts1Cje Mouse Model. *Am J Hum Genet.* doi: 10.1016/j.ajhg.2020.10.001.
29. Moustafa, A. et al. (2020). Changes in nitric oxide, carbon monoxide, hydrogen sulfide and male reproductive hormones in response to chronic restraint stress in rats. *Free Radic Biol Med.* doi: 10.1016/j.freeradbiomed.2020.10.315.
30. Palani, A. et al (2020). Impact of oxidative stress on semen parameters in normozoospermic infertile men: a case-control study. *Afr J Urol.* doi: 10.1186/s12301-020-00061-6.

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's 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

©2010-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.