

---

Product Manual

# OxiSelect™ TBARS Assay Kit (MDA Quantitation), Trial Size

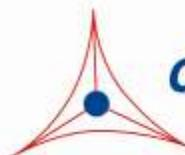
Catalog Number

STA-330-T

20 assays

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

---



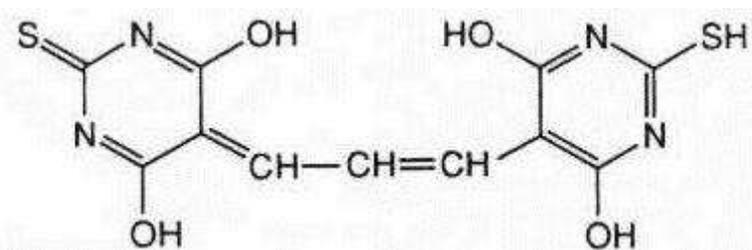
**CELL BIOLABS, INC.**

*Creating Solutions for Life Science Research*

## **Introduction**

Lipid peroxidation is a well-defined mechanism of cellular damage in animals and plants. Lipid peroxides are unstable indicators of oxidative stress in cells that decompose to form more complex and reactive compounds such as Malondialdehyde (MDA) and 4-hydroxynonenal (4-HNE), natural by-products of lipid peroxidation. Oxidative modification of lipids can be induced *in vitro* by a wide array of pro-oxidant agents and occurs *in vivo* during aging and in certain disease conditions. Measuring the end products of lipid peroxidation is one of the most widely accepted assays for oxidative damage. These aldehydic secondary products of lipid peroxidation are generally accepted markers of oxidative stress.

Thiobarbituric Acid Reactive Substances (TBARS) is a well-established assay for screening and monitoring lipid peroxidation. The rapid and easy protocol has been modified by researchers in the evaluation of drugs, food, as well as human and animal tissue samples. MDA forms a 1:2 adduct with thiobarbituric acid (Figure 1). The MDA-TBA adduct formed from the reaction of MDA in samples with TBA can be measured colorimetrically or fluorometrically. TBARS levels are determined from a Malondialdehyde equivalence standard.



**Figure 1. MDA-TBA Adduct**

The TBARS Assay has provided relevant information concerning free radical activity in disease states and measurement of many compounds anti-oxidant characteristics. Although the specificity of TBARS toward compounds other than MDA has been controversial, the assay continues to be the most widely employed format for monitoring lipid peroxidation. Lipids with higher degrees of unsaturated bonds produce higher TBARS values. Interfering soluble TBARS can be minimized if lipoprotein fractions are first acid precipitated from samples. Biological samples may contain a mixture of thiobarbituric acid reactive substances such as hydroperoxides and aldehydes, which increase in response to oxidative stress. If excessively high TBARS values are obtained, a more specific assay such as HPLC should be employed.

The OxiSelect™ TBARS Assay Kit offers a simple, reproducible, and consistent system for the detection of lipid peroxidation in urine, plasma, serum, lysates, and tissue homogenates. This kit includes an MDA standard for use as a positive control. Each Trial Size TBARS Assay Kit provides sufficient reagents to perform 20 tests including standard curve and unknown samples.

## **Assay Principle**

The Thiobarbituric Acid Reactive Substances (TBARS) Assay Kit is a tool for the direct quantitative measurement of MDA in biological samples. The unknown MDA containing samples or MDA standards are first reacted with TBA at 95°C. After a brief incubation, the samples and standards can

be read either spectrophotometrically or fluorometrically. The MDA content in unknown samples is determined by comparison with the predetermined MDA standard curve.

### **Related Products**

1. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
2. STA-325: OxiSelect™ Oxidative RNA Damage ELISA Kit (8-OHG Quantitation)
3. STA-333: MDA-BSA Control
4. STA-337: OxiSelect™ 8-iso-Prostaglandin F2a Activity Assay Kit
5. STA-344: OxiSelect™ Hydrogen Peroxide/Peroxidase Assay Kit
6. STA-347: OxiSelect™ In Vitro ROS/RNS Assay Kit (Green Fluorescence)
7. STA-816: OxiSelect™ N-epsilon-(Carboxymethyl) Lysine (CML) Competitive ELISA Kit
8. STA-817: OxiSelect™ Advanced Glycation End Products (AGE) Competitive ELISA Kit
9. STA-832: OxiSelect™ MDA Adduct Competitive ELISA Kit
10. STA-838: OxiSelect™ HNE Adduct Competitive ELISA Kit

### **Kit Components**

1. MDA Standard (Part No. 233001-T): One 125 µL amber vial of 1.0 mM Malondialdehyde bis(dimethyl acetal)
2. Thiobarbituric Acid (TBA) (Part No. 233002-T): One 100 mg tube
3. SDS Lysis Solution (Part No. 233003-T): One 2 mL tube
4. 2X TBA Acid Diluent (Part No. 233004-T): Two 1.5 mL tubes
5. Sodium Hydroxide Solution (Part No. 233005-T): One 0.5 mL tube
6. 100X BHT Solution (Part No. 233006-T): One 100 µL tube of 5% Butylated hydroxytoluene (BHT) in methanol

### **Materials Not Supplied**

1. MDA samples: plasma, serum, urine, tissue or cell lysate
2. 1X PBS
3. Microcentrifuge and conical tubes
4. Heat block, incubator or water bath
5. n-Butanol
6. 96-well clear, flat-bottomed microplate for reading samples/standards
7. 96-well black fluorescence microplate for reading samples/standards
8. Spectrophotometric microplate reader capable of reading at 532nm
9. Fluorometric microplate reader capable of reading at 540nm excitation and 590nm emission

## **Storage**

Store all components at 4°C.

## **Preparation of Reagents**

- 1X TBA Acid Diluent: Dilute the 2X TBA Acid Diluent with equal parts distilled or deionized water.
- SDS Lysis Solution: If precipitated crystals are present, briefly heat the solution at 37°C to redissolve the SDS crystals.
- TBA Reagent: Prepare the TBA Reagent just before use. Prepare a 5.2 mg/mL solution of TBA Reagent by weighing out an amount of TBA needed for all samples and standards (eg: 13 mg of TBA is enough to prepare 10 tests). Add 1X TBA Acid Diluent (see above) to the TBA and stir or mix vigorously until the powder has dissolved (e.g: 2.5 mL 1X TBA Diluent for 13 mg of TBA). Adjust the pH of the solution to pH 3.5 with the Sodium Hydroxide Solution.

*Note: The TBA Reagent is stable for 24 hours. Do not store or reuse diluted solutions.*

## **Preparation of Samples**

***Important Note: All samples should be assayed immediately upon collection or stored at -80°C for up to 1-2 months.***

- Tissue: Because hemoglobin interferes with the assay, blood should be removed from tissue sample by perfusion with PBS containing heparin. Resuspend tissue at 50 to 100 mg/mL in PBS. To prevent further oxidation, add 100X BHT Solution to achieve a final concentration of 1X (for example, add 10 µL of 100X BHT to 1 mL of sample volume). Homogenize the tissue sample on ice, spin at 10,000 g for 5 min to collect the supernatant. The supernatant can be assayed directly for its TBARS level and results can be normalized based on its protein concentration.
- Plasma: To minimize the hemoglobin interference, prepare the plasma sample as soon as possible after blood being drawn. To prevent further oxidation, add 100X BHT to plasma samples to achieve a final concentration of 1X (for example, add 10 µL of 100X BHT to 1 mL of plasma). Plasma samples can be assayed directly without further processing.
- Cells: Resuspend cells at 1-2 x 10<sup>7</sup> cells/mL in PBS. To prevent further oxidation, add 100X BHT Solution to achieve a final concentration of 1X (for example, add 10 µL of 100X BHT to 1 mL of cell suspension). Homogenize or sonicate the cells on ice. Use the whole homogenate in the assay.
- Urine: To remove insoluble particles, spin at 10,000 g for 5 min. The supernatant can be assayed directly.

## **Preparation of Standard Curve**

Prepare a dilution series of MDA standards in the concentration range of 125  $\mu\text{M}$  – 0  $\mu\text{M}$  by diluting the MDA Standard in distilled or deionized water (Table 1). It is recommended that standards be performed in duplicate.

<b>Standard Tubes</b>	<b>MDA Standard (<math>\mu\text{L}</math>)</b>	<b>Water (<math>\mu\text{L}</math>)</b>	<b>MDA Standard (<math>\mu\text{M}</math>)</b>
1	75 $\mu\text{L}$	525 $\mu\text{L}$	125
2	250 $\mu\text{L}$ of Tube #1	250 $\mu\text{L}$	62.5
3	250 $\mu\text{L}$ of Tube #2	250 $\mu\text{L}$	31.25
4	250 $\mu\text{L}$ of Tube #3	250 $\mu\text{L}$	15.63
5	250 $\mu\text{L}$ of Tube #4	250 $\mu\text{L}$	7.81
6	250 $\mu\text{L}$ of Tube #5	250 $\mu\text{L}$	3.91
7	250 $\mu\text{L}$ of Tube #6	250 $\mu\text{L}$	1.95
8	250 $\mu\text{L}$ of Tube #7	250 $\mu\text{L}$	0.98
9	0 $\mu\text{L}$	250 $\mu\text{L}$	0.0

**Table 1. Preparation of MDA Standards**

## **Assay Protocol**

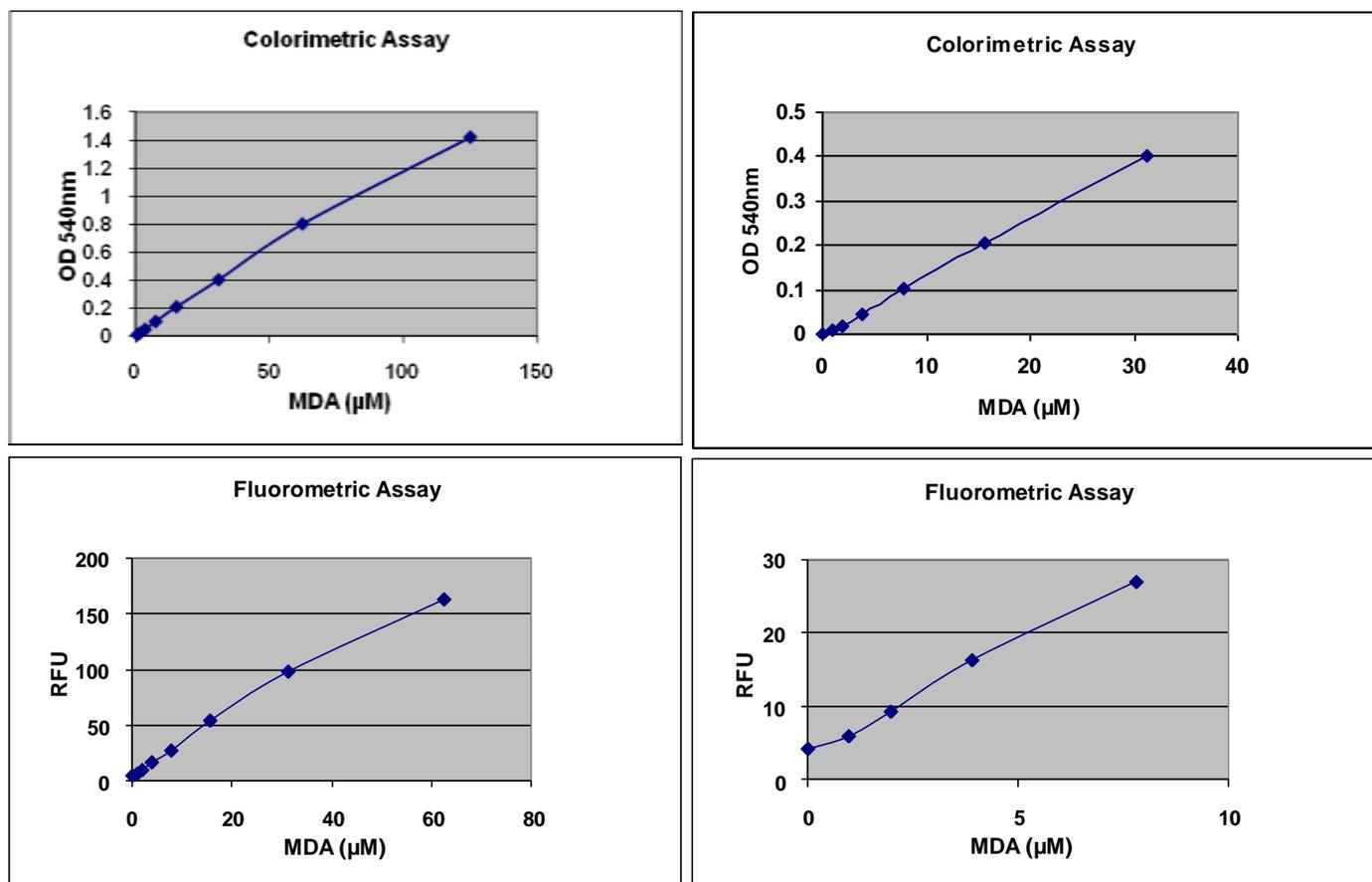
1. Prepare and mix all reagents thoroughly before use. Each MDA-containing sample and standard should be assayed in duplicate. High content MDA samples can be further diluted for analysis.
2. Add 100  $\mu\text{L}$  of unknown samples or MDA standards to separate microcentrifuge tubes.
3. Add 100  $\mu\text{L}$  of the SDS Lysis Solution to both the unknown samples and the MDA standards. Mix thoroughly. Incubate samples for 5 minutes at room temperature.
4. Add 250  $\mu\text{L}$  of TBA Reagent to each sample and standard to be tested.
5. Close each tube and incubate at 95°C for 45-60 minutes.
6. Remove tubes and cool to room temperature in an ice bath for 5 minutes.
7. Centrifuge all sample tubes at 3000 rpm for 15 minutes. Remove the supernatant from samples for further analysis.
8. (optional) Butanol Extraction: To prevent the interference of hemoglobin and its derivatives, we recommend the following extraction procedure:
  - a. Transfer 300  $\mu\text{L}$  of the supernatant (Step 7) to another tube, add 300  $\mu\text{L}$  of n-Butanol. Vortex vigorously for 1-2 minutes and centrifuge for 5 minutes at 10,000 g.
  - b. Transfer the butanol fraction for further measurement.
9. Spectrophotometric Measurement: Transfer 200  $\mu\text{L}$  of the MDA standards and samples to a 96 well microplate compatible with a spectrophotometric plate reader. Remember to include a 0  $\mu\text{M}$

blank control. It is recommended that duplicates of each standard and sample should be read. Read the absorbance at 532nm.

**Fluorometric Measurement:** Transfer 150  $\mu\text{L}$  of the MDA standards and samples to a 96 well black fluorescence microplate compatible with a fluorometric plate reader. Remember to include a 0  $\mu\text{M}$  blank control. It is recommended that duplicates of each standard and sample should be read. Read the plate at 540 nm excitation and 590 nm emission.

### **Example of Results**

The following figures demonstrate typical MDA Quantitation results by TBARS Assay Kit. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 2. MDA Standard Curve.** The MDA standard curve was created as described in the Assay Protocol. Top Panel: Colorimetric Detection; Bottom Panel: Fluorometric Detection.

### **References**

1. Armstrong, D. and Browne, R. (1994). *Free Radicals in Diagnostic Medicine*. 366: 43-58.
2. Armstrong, D., et al. (1998). *Free Radicals and Antioxidant Protocols*. 108: 315-324.
3. Boyum, A. (1966). *J. of Clinical Investigation*. 21: Supplement 97.
4. Braun, D. and Fromherz, P. (1997). *Applied Physics A*.

5. Gidez, L., et al. (1982). *J. of Lipid Research*. 23: 1206-1223.
6. Lef'evre G., et al. (1998). *Annals de Biologie Clinique*. 56(3): 305-319.
7. Ohkawa, H., et al. (1979). *Anal. Biochem*. 95: 351-358.
8. Yagi, K. (1998). *Free Radicals and Antioxidant Protocols*. 108: 101-106.

### **Recent Product Citations**

1. 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.
2. Bhattarai, K.R. et al. (2021). TMBIM6 regulates redox-associated posttranslational modifications of IRE1 $\alpha$  and ER stress response failure in aging mice and humans. *Redox Biol.* **47**:102128. doi: 10.1016/j.redox.2021.102128.
3. Zhao, X. & Chen, F. et al. (2021). Propofol induces the ferroptosis of colorectal cancer cells by downregulating STAT3 expression. *Oncol Lett.* **22**(5):767. doi: 10.3892/ol.2021.13028.
4. Raj, P. et al. (2021). Saskatoon berry supplementation prevents cardiac remodeling without improving renal disease in an animal model of reno-cardiac syndrome. *J Food Biochem*. doi: 10.1111/jfbc.13893.
5. Liang, Y. et al. (2021). Oxidative stress and EROD activity in Caco-2 cells upon exposure to chlorinated hydrophobic organic compounds from drinking water reservoirs. *Sci Total Environ.* **804**:150015. doi: 10.1016/j.scitotenv.2021.150015.
6. Aldayel, T.S. et al. (2021). Assessment of modulatory activity of *Uncaria tomentosa* extract against fipronil immunotoxicity in male rats. *Ecotoxicol Environ Saf.* **224**:112674. doi: 10.1016/j.ecoenv.2021.112674.
7. Ammar, I.M.M. & Salem, M.A.A. et al. (2021). Amelioration of polycystic ovary syndrome-related disorders by supplementation of thymoquinone and metformin. *Middle East Fertil Soc J.* doi: 10.1186/s43043-021-00076-1.
8. Martínez, M.A. et al. (2021). Brown marine algae *Gongolaria baccata* extract protects Caco-2 cells from oxidative stress induced by tert-butyl hydroperoxide. *Food Chem Toxicol.* doi: 10.1016/j.fct.2021.112460.
9. 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.anifeedsci.2021.114994.
10. 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.
11. Kim, J.U. et al. (2021). Inactivation of *Bacillus cereus* spores using a combined treatment of UV-TiO<sub>2</sub> photocatalysis and high hydrostatic pressure. *Innov Food Sci Emerg Technol.* doi: 10.1016/j.ifset.2021.102676.
12. 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.bbali.2021.158927.
13. Zhou, Z. et al. (2021). Exogenous activation of toll-like receptor 5 signaling mitigates acetaminophen-induced hepatotoxicity in mice. *Toxicol Lett.* doi: 10.1016/j.toxlet.2021.01.022.

14. Yu, C. et al. (2021). Ferulic acid exerts Nrf2-dependent protection against prenatal lead exposure-induced cognitive impairment in offspring mice: Running title: Cognition-enhancing effects of ferulic acid. *J Nutr Biochem*. doi: 10.1016/j.jnutbio.2021.108603.
15. McCollum, C.R. et al. (2021). Photoexcited Quantum Dots as Efficacious and Nontoxic Antibiotics in an Animal Model. *ACS Biomater Sci Eng*. doi: 10.1021/acsbiomaterials.0c01406.
16. Huang, Y. et al. (2020). Cataract formation in transgenic HO-1 G143H mutant mice: Involvement of oxidative stress and endoplasmic reticulum stress. *Biochem Biophys Res Commun*. **537**:43-49. doi: 10.1016/j.bbrc.2020.12.071.
17. 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.
18. Wang, J. et al. (2021). Cardiorespiratory responses in healthy young adults with exposure to indoor airborne PAEs: A randomized, crossover trial of air purification. *Environ Int*. **156**:106761. doi: 10.1016/j.envint.2021.106761.
19. Yuan, X. et al. (2021). Toxicometabolomics of lindane in adult zebrafish (*Danio rerio*) using GC-MS/MS and LC-Orbitrap-MS/MS. *Appl Biol Chem*. doi: 10.1186/s13765-021-00623-4.
20. Lee, J.W. et al. (2021). Protective Effects of Topical Application of Nitrite on Testicular Ischemia-Reperfusion Injury in Rats. *Oxid Med Cell Longev*. doi: 10.1155/2021/5514537.
21. Toriumi, K. et al. (2021). Combined glyoxalase 1 dysfunction and vitamin B6 deficiency in a schizophrenia model system causes mitochondrial dysfunction in the prefrontal cortex. *Redox Biol*. **45**:102057. doi: 10.1016/j.redox.2021.102057.
22. Duarte, M.E. et al. (2021). Modulation of jejunal mucosa-associated microbiota in relation to intestinal health and nutrient digestibility in pigs by supplementation of  $\beta$ -glucanase to corn soybean meal-based diets with xylanase. *J Anim Sci*. doi: 10.1093/jas/skab190.
23. Shen, Y.B. et al. (2021). Physiological Effects of Deoxynivalenol from Naturally Contaminated Corn on Cerebral Tryptophan Metabolism, Behavioral Response, Gastrointestinal Immune Status and Health in Pigs Following a Pair-Feeding Model. *Toxins*. **13**(6):393. doi: 10.3390/toxins13060393.
24. Lee, S.H. et al. (2021). Dietary soluble flaxseed oils as a source of omega-3 polyunsaturated fatty acids for laying hens. *Poult Sci*. doi: 10.1016/j.psj.2021.101276.
25. Alsas, M. & Alzeer, S. (2021). Evaluation of some biological parameters of gasoline station attendants in Damascus, Syria. *Heliyon*. **7**(5):e07056. doi: 10.1016/j.heliyon.2021.e07056.
26. Malicka, M. et al. (2021). Differences in the effects of single and mixed species of AMF on the growth and oxidative stress defense in *Lolium perenne* exposed to hydrocarbons. *Ecotoxicol Environ Saf*. **217**:112252. doi: 10.1016/j.ecoenv.2021.112252.
27. Malcok, U.A. et al. (2021). Protective effect of metformin against lithium-induced cerebral neurotoxicity in rats. *Med Science*. **10**(2):350-5. doi: 10.5455/medscience.2020.12.253.
28. Lauridsen, C. et al. (2021). Effect of Maternal Dietary Redox Levels on Antioxidative Status and Immunity of the Suckling Off-Spring. *Antioxidants (Basel)*. **10**(3):478. doi: 10.3390/antiox10030478.
29. Sun, Y. et al. (2021). Dietary inclusion of multispecies probiotics to reduce the severity of post-weaning diarrhea caused by F18+ *E. coli* in pigs. *Anim Nutr*. doi: 10.1016/j.aninu.2020.08.012.
30. Sangweni, N.F. et al. (2021). The triterpene, methyl-3 $\beta$ -hydroxylanosta-9,24-dien-21-oate (RA3), attenuates high glucose-induced oxidative damage and apoptosis by improving energy metabolism. *Phytomedicine*. doi: 10.1016/j.phymed.2021.153546.

## **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](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2013-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.