

---

Product Manual

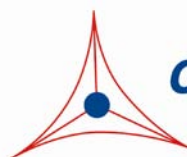
# OxiSelect™ Protein Carbonyl ELISA Kit

## Catalog Number

STA-310	96 assays
STA-310-5	5 x 96 assays

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

---



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

## **Introduction**

Protein oxidation is defined as the covalent modification of a protein induced either directly by reactive oxygen species (ROS) or indirectly by reaction with secondary by-products of oxidative stress. Oxidative modification of proteins can be induced in vitro by a wide array of pro-oxidant agents and occurs in vivo during aging and in certain disease conditions.

There are numerous types of protein oxidative modification. The most common products of protein oxidation in biological samples are the protein carbonyl derivatives of Pro, Arg, Lys, and Thr. These derivatives are chemically stable and serve as markers of oxidative stress for most types of ROS. Many of the current assays involve derivatization of the carbonyl group with dinitrophenylhydrazine (DNPH), followed by immunoblotting with an anti-DNP antibody. The Protein Carbonyl ELISA was first developed by Buss and co-workers, the protein samples ( $\geq 4$  mg/mL) react with DNPH and then adsorb to wells of an ELISA plate before probe with anti-DNPH antibody. In their method, protein samples containing low amounts of protein must be concentrated to at least 4 mg/mL by TCA precipitation. However, TCA precipitation results a 20% loss of the total carbonyl values, and loss of protein during precipitation is also expected. In Cell Biolabs' OxiSelect™ Protein Carbonyl ELISA Kit, protein samples are first allowed to adsorb to wells of a 96-well plate and then react with DNPH. There is no need to concentrate protein in experimental and clinical samples with low amounts of protein ( $< 4$  mg/mL) and the kit requires protein sample as little as 10  $\mu$ g/mL.

The OxiSelect™ Protein Carbonyl ELISA Kit is an enzyme immunoassay developed for rapid detection and quantitation of protein carbonyls. The quantity of protein carbonyls in protein sample is determined by comparing its absorbance with that of a known reduced/oxidized BSA standard curve. Each kit provides sufficient reagents to perform up to 96 assays, including standard curve and unknown protein samples.

## **Assay Principle**

BSA standards or protein samples (10  $\mu$ g/mL) are adsorbed onto a 96-well plate for 2 hrs at 37°C. The protein carbonyls present in the sample or standard are derivatized to DNP hydrazone and probed with an anti-DNP antibody, followed by an HRP conjugated secondary antibody. The protein carbonyl content in unknown sample is determined by comparing with a standard curve that is prepared from predetermined reduced and oxidized BSA standards.

## **Related Products**

1. STA-303: OxiSelect™ Nitrotyrosine Protein Immunoblot Kit
2. STA-304: Protein Tyrosine Nitration Control (Nitrotyrosine-BSA)
3. STA-305: OxiSelect™ Nitrotyrosine ELISA Kit
4. STA-308: OxiSelect™ Protein Carbonyl Immunoblot Kit
5. STA-309: Oxidized Protein Immunoblot Control (Carbonyl-BSA)
6. STA-318: OxiSelect™ AOPP Assay Kit
7. STA-319: AOPP-Human Serum Albumin (AOPP-HSA)

8. STA-816: OxiSelect™ N-epsilon-(Carboxymethyl) Lysine (CML) Competitive ELISA Kit
9. STA-817: OxiSelect™ Advanced Glycation End Products (AGE) Competitive ELISA Kit

### **Kit Components**

1. 96-well Protein Binding Plate (Part No. 231001): One strip well 96-well plate.
2. Anti-DNP Antibody (1000X) (Part No. 231002): One 20 µL vial of anti-DNP Rabbit IgG.
3. Secondary Antibody, HRP Conjugate (1000X) (Part No. 231009): One 20 µL vial.
4. DNPH (Part No. 231004): One 200 mg amber vial.
5. 2X DNPH Diluent (Part No. 231005): One 15 mL bottle.
6. Blocking Reagent (Part No. 231006): One 20 g bottle.
7. 10X Wash Buffer (Part No. 310806): One 100 mL bottle.
8. Substrate Solution (Part No. 310807): One 12 mL amber bottle.
9. Stop Solution (Part. No. 310808): One 12 mL bottle.
10. Reduced BSA Standard (Part No. 231007): One 200 µL vial of 1 mg/mL fully reduced BSA in PBS.
11. Oxidized BSA Standard (Part No. 231008): One 200 µL vial of 1 mg/mL oxidized BSA in PBS at 7.5 nmol protein carbonyl/mg proteins. The protein carbonyl is predetermined by a spectrophotometric method as described by Reznick and Parker (See Ref. 5).

### **Materials Not Supplied**

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

### **Storage**

Upon receipt, aliquot and store both the Reduced and Oxidized BSA Standards at -20°C to avoid multiple freeze/thaw cycles. Store all other components at 4°C.

### **Preparation of Reagents**

- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.
- 1X DNPH Diluent: Dilute the 2X DNPH Diluent to 1X with deionized water. Mix well.

- Blocking Solution: Weigh out 5 g of Blocking Reagent, dissolve in 100 mL of 1X PBS, and store at 4°C for up to one week.
- DNPH Working Solution: Weigh out 1-5 mg of DNPH, and dissolve in 1X DNPH diluent to 1 mg/mL. This DNPH stock solution is stable for one week when stored in the dark at 4°C. Based on the number of tests, FRESHLY prepare appropriate amount of DNPH Working Solution by diluting the 1 mg/mL stock DNPH solution to 0.04 mg/mL in 1X DNPH diluent. Example: for 20 assays, transfer 100  $\mu$ L of 1 mg/mL DNPH stock solution to a tube containing 2.4 mL of 1X DNPH Diluent, mix well and use it IMMEDIATELY.
- Anti-DNP Antibody and Secondary Antibody: Immediately before use dilute the Anti-DNPH antibody 1:1000 and Secondary Antibody 1:1000 with 1X Blocking Solution. Do not store diluted solutions.

### **Preparation of Protein Carbonyl BSA Standards**

1. Freshly Prepare 10  $\mu$ g/mL of reduced or oxidized BSA by diluting the 1 mg/mL BSA standards in 1X PBS. Example: Add 20  $\mu$ L to 1.98 mL of 1X PBS.
2. Prepare a series of carbonyl BSA standards by mixing the oxidized BSA and reduced BSA in the proper ratios according to Table 1.

Standard Tubes	10 $\mu$ g/mL Oxidized BSA ( $\mu$ L)	10 $\mu$ g/mL Reduced BSA ( $\mu$ L)	[Protein Carbonyl] (nmol/mg)
1	400	0	7.5
2	320	80	6.0
3	240	160	4.5
4	160	240	3.0
5	80	320	1.5
6	40	360	0.75
7	20	380	0.375
8	0	400	0

**Table 1. Preparation of Protein Carbonyl BSA Standard Curve**

### **Preparation of Samples**

1. Perform a protein assay such as Bradford or BCA on all samples to determine the protein concentration.

*Notes for cell and tissue lysates:*

- Lysates should not be prepared in lysis buffer containing Triton X-100, NP-40, or Igepal CA-630 because these detergents interfere with protein coating of the plate unless the detergent concentration in the 10  $\mu$ g/mL protein samples is no more than 0.001%. We recommend lysis by homogenization or sonication.
- A high concentration of nucleic acid in cell or tissue lysates can erroneously contribute to higher estimation of carbonyl content. To remove nucleic acid, we recommend one of the following procedures:

- 1. Pretreat lysate with nuclease, followed by ammonium sulfate precipitation of high percentage saturation.*
- 2. Add streptomycin sulfate or PEI to a final concentration of 1% and 0.5% respectively, incubate 30 minutes at room temperature and remove the nuclei acid precipitates by centrifuging at 6000 g for 10 minutes at 4°C.*

2. Dilute each protein sample to 10 µg/mL in 1X PBS prior to use in the assay.

*Note: Samples with high concentrations of protein carbonyl content may be further diluted 5-10 fold in 10 µg/mL Reduced BSA. A titration may be performed to ensure the samples fall in the range of the standard curve.*

### **Assay Protocol**

1. Prepare unknown samples according to the Preparation of Samples section above. Each 10 µg/mL protein sample and BSA Standard should be assayed in duplicate or triplicate.
2. Add 100 µL of 10 µg/mL protein samples or reduced/oxidized BSA standards to the 96-well Protein Binding Plate. Incubate at 37°C for at least 2 hours or 4°C overnight.
3. Wash wells 3 times with 250 µL 1X PBS per well. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess wash solution.
4. Add 100 µL of the DNPH Working Solution and incubate for 45 minutes at room temperature in the dark.
5. Wash wells with 250 µL of 1X PBS/Ethanol (1:1, v/v) with incubation on an orbital shaker for 5 minutes. Repeat washing a total of 5 times, aspirating between each. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess wash solution. Wash 2 times with 250 µL of 1X PBS.
6. Add 200 µL of Blocking Solution per well and incubate for 1-2 hours at room temperature on an orbital shaker.
7. Wash 3 times with 250 µL of 1X Wash Buffer 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.
8. Add 100 µL of the diluted anti-DNP antibody to all wells and incubate for 1 hour at room temperature on an orbital shaker. Wash the strip wells 3 times according to step 7 above.
9. Add 100 µL of the diluted HRP conjugated secondary antibody to all wells and incubate for 1 hour at room temperature on an orbital shaker. Wash the strip wells 5 times according to step 7 above.

10. Warm Substrate Solution to room temperature. Add 100  $\mu$ L 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.

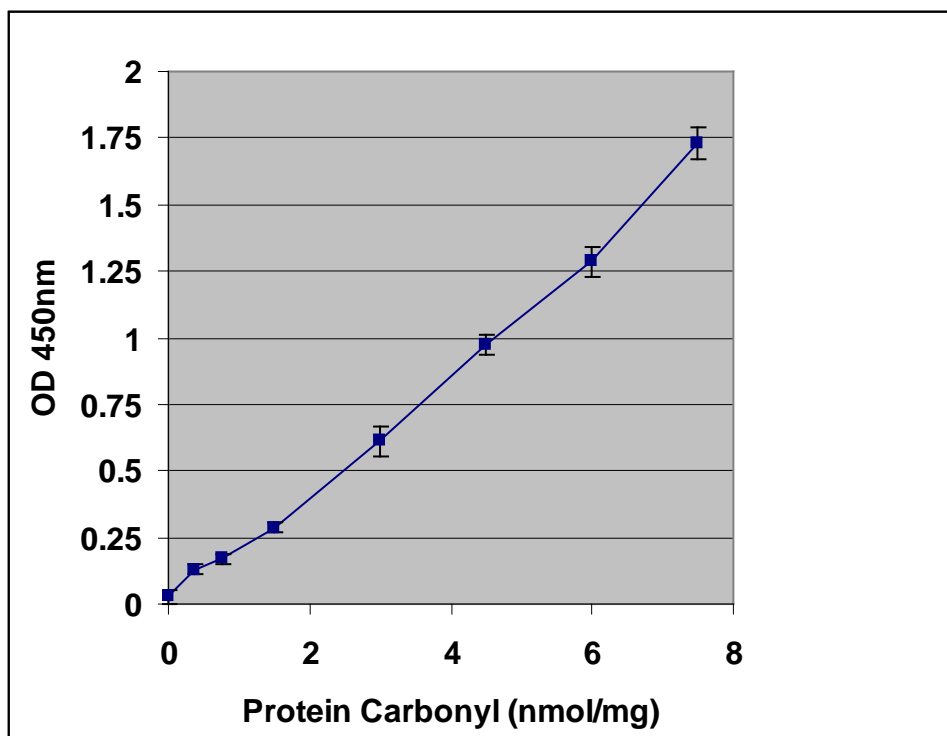
*Note: Watch plate carefully; if color changes rapidly, the reaction may need to be stopped sooner to prevent saturation.*

11. Stop the enzyme reaction by adding 100  $\mu$ L of Stop Solution to each well. Results should be read immediately (color will fade over time).

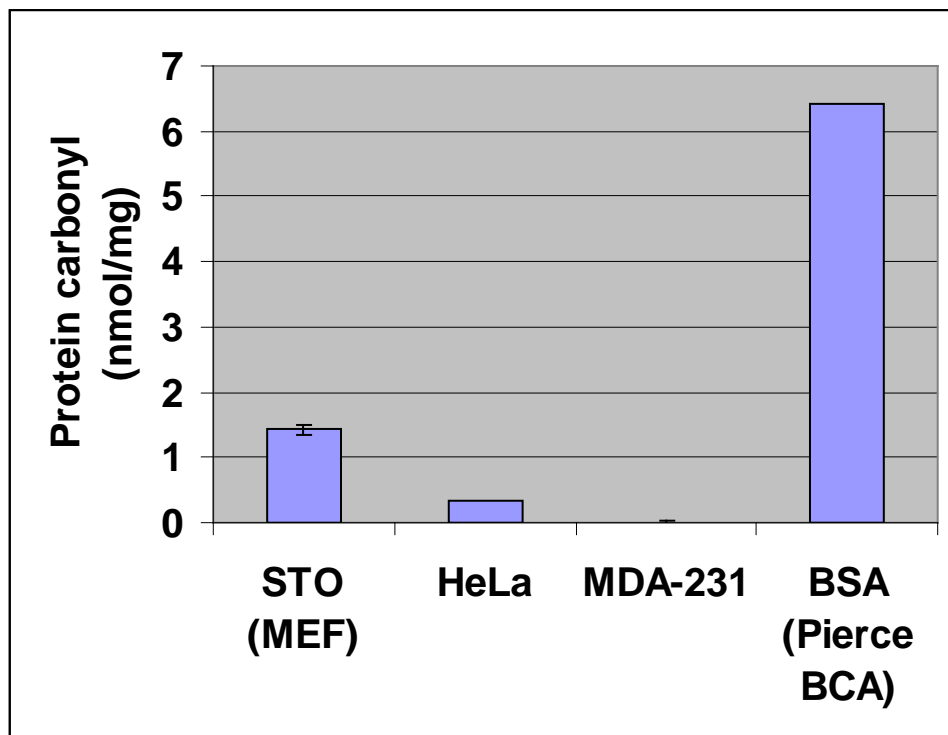
12. Read absorbance of each well on a plate reader using 450 nm as the primary wave length. Using the fully reduced BSA standard as absorbance blank.

### **Example of Results**

The following figures demonstrate typical Protein Carbonyl ELISA results. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1: Protein Carbonyl ELISA Standard Curve**



**Figure 2: Amount of Protein Carbonyl Content for Cell Lysate and BSA Standard.** STO (MEF), HeLa and MDA-231 cells were sonicated in 25mM HEPES, pH 7.5, 150 mM NaCl, 10 mM MgCl<sub>2</sub>, 1 mM EDTA, 2% Glycerol. Cell Lysates and BSA Standard from Pierce BCA Protein Assay were diluted to 10 µg/mL with 1X PBS and coated onto a 96-well Protein Binding Plate. The protein carbonyl levels were determined as described in the Assay Protocol.

## References

1. Cadenas, E., Boveris, A., Ragan, CI., and Stoppani, AO. (1977). *Archives of Biochemistry & Biophysics* **180**:248–257.
2. Wakeyama, H., Takeshige, K., Takayanagi, R., and Minakami, S. (1982). *Biochem J.* **205**:593–601.
3. Talent, JM., Kong, Y., and Gracy, RW. (1998). *Anal. Biochem.* **263**:31–38.
4. Buss, H, Chan, TP, Sluis, KB, Domigan, NM, and Winterbourn, CC. (1985) *Free Radic Biol Med.* **23**:361-6.
5. Reznick, AZ., and Packer, L. (1994) *Methods Enzymol.* **233**:263-357.

## Recent Product Citations

1. Ren, Z., et al. (2017). Ethanol-induced Damage to the Developing Spinal Cord: The Involvement of CCR2 Signaling. *Biochim Biophys Acta*. pii: S0925-4439(17)30270-3. doi: 10.1016/j.bbadis.2017.07.035.
2. Hondur G, et al. (2017). Oxidative Stress-Related Molecular Biomarker Candidates for Glaucoma. *Invest Ophthalmol Vis Sci.* 58(10):4078-4088. doi: 10.1167/iovs.17-22242.
3. Kim, L., et al. (2017). Antioxidant activities of ethanolic and acidic ethanolic extracts of astringent persimmon in H<sub>2</sub>O<sub>2</sub>-stimulated Caco-2 human colonic epithelial cells. *Food Sci Biotechnol.* **26**: 1085.

4. Fan, N. et al. (2017). Rapid repeatable in vivo detection of retinal reactive oxygen species. *Exp. Eye Res.* **161**:71-81.
5. Strickland, F.M. et al. (2017). Oxidative T cell modifications in lupus and Sjogren's syndrome. *Lupus* **2(1)**:121.
6. Fortin, E. et al. (2017). Linoleic acid,  $\alpha$ -linolenic acid and enterolactone affect lipid oxidation and expression of lipid metabolism and antioxidant-related genes in hepatic tissue of dairy cows. *Br. J. Nutr.* **117(9)**:1199-1211.
7. Kawai, T. et al. (2017). Unconventional role of voltage-gated proton channels (VSOP/Hv1) in regulation of microglial ROS production. *J. Neurochem.* doi:10.1111/jnc.14106.
8. Tangtrongsup, S. and Kisiday, J.D. (2017). Modulating the oxidative environment during mesenchymal stem cells chondrogenesis with serum increases collagen accumulation in agarose culture. *J. Orthop. Res.* doi:10.1002/jor.23618.
9. Bulteau A.L., et al. (2017). Dysfunction of mitochondrial Lon protease and identification of oxidized protein in mouse brain following exposure to MPTP: Implications for Parkinson disease. *Free Radic Biol Med.* **108**:236-246. doi: 10.1016/j.freeradbiomed.2017.03.036.
10. Zhang, M. et al. (2017). Allicin Decreases Lipopolysaccharide-Induced Oxidative Stress and Inflammation in Human Umbilical Vein Endothelial Cells through Suppression of Mitochondrial Dysfunction and Activation of Nrf2. *Cell Physiol. Biochem.* **41(6)**:2255-2267.
11. Ahn, M.Y., et al. (2017). Anti-aging effect and gene expression profiling of dung beetle glycosaminoglycan in aged rats. *Biomater Res.* **21**:5. doi: 10.1186/s40824-017-0091-9.
12. Lundh, M., et al. (2017). The immunoproteasome is induced by cytokines and regulates apoptosis in human islets. *J Endocrinol.* pii: JOE-17-0110. doi: 10.1530/JOE-17-0110.
13. Bulteau, A.L., et al. (2017). Dysfunction of mitochondrial Lon protease and identification of oxidized protein in mouse brain following exposure to MPTP: Implications for Parkinson disease. *Free Radic Biol Med.* **108**:236-246. doi: 10.1016/j.freeradbiomed.2017.03.036.
14. Voutetakis, K. et al. (2017). Anti-ageing properties of Khelma Longevity™: treatment of human fibroblasts increases proteasome levels and decreases the levels of oxidized proteins. *N Biotechnol.* pii: S1871-6784(17)30134-6. doi: 10.1016/j.nbt.2017.03.002.
15. Francioli, C. et al. (2017). Pyrrolidine dithiocarbamate administered during ex-vivo lung perfusion promotes rehabilitation of injured donor rat lungs obtained after prolonged warm ischemia. *PLoS One.* **12(3)**:e0173916. doi: 10.1371/journal.pone.0173916.
16. Ahn, M.Y. et al. (2017). Anti-Obesity Effect of *Bombus ignitus* Queen Glycosaminoglycans in Rats on a High-Fat Diet. *Int J Mol Sci.* **18(3)**. pii: E681. doi: 10.3390/ijms18030681.
17. Kim, B.G. et al. (2017). Effect of TiO<sub>2</sub> Nanoparticles on Inflammation-Mediated Airway Inflammation and Responsiveness. *Allergy Asthma Immunol Res.*
18. Kim, E.C. et al (2017). Screening and Characterization of Drugs That Protect Corneal Endothelial Cells Against Unfolded Protein Response and Oxidative Stress. *Invest. Ophthalmol. Vis. Sci.* **58(2)**:892-900. doi: 10.1167/iovs.16-20147.
19. Mostek, A. et al. (2017). Cryopreservation of bull semen is associated with carbonylation of sperm proteins. *Theriogenology.* **92**:95-102. doi: 10.1016/j.theriogenology.2017.01.011.
20. Stadem, P.S. et al. (2017). Markers of oxidative stress in umbilical cord blood from G6PD deficient African newborns. *PLoS One.* **12(2)**:e0172980. doi: 10.1371/journal.pone.0172980.
21. Wang, N. et al. (2017). Apigenin attenuates streptozotocin-induced pancreatic  $\beta$  cell damage by its protective effects on cellular antioxidant defense. *In Vitro Cell.Dev.Biol.Animal.* doi:10.1007/s11626-017-0135-4.



22. Webb, S.J. et al. (2016). Stress physiology and weapon integrity of intertidal mantis shrimp under future ocean conditions. *Sci. Rep.* **6**:38637.
23. Xia, M. et al. (2016). Remote limb ischemic preconditioning protects rats against cerebral ischemia via HIF-1 $\alpha$ /AMPK/HSP70 pathway. *Cell. Mol. Neurobiol.* doi:10.1007/s10571-016-0444-2.
24. Augustyniak, E. et al. (2016). Validation of protein carbonyl measurement: a multi-centre study. *Redox Biol.* **4**:149-157.
25. Sakellariou, G. K. et al. (2016). Mitochondrial ROS regulate oxidative damage and mitophagy but not age-related muscle fiber atrophy. *Sci. Rep.* doi:10.1038/srep33944.
26. Choi, Y. S. et al. (2016). Preservation of myocardial fatty acid oxidation prevents diastolic dysfunction in mice subjected to angiotensin II infusion. *J Mol Cell Cardiol.* doi:10.1016/j.yjmcc.2016.09.001.
27. Li, Y. et al. (2016). Visible-light-driven photocatalytic inactivation of MS2 by metal-free gC 3 N 4: Virucidal performance and mechanism. *Water Res.* doi:10.1016/j.watres.2016.10.009.
28. Laubertova, L. et al. (2016). Oxidative stress markers and their dynamic changes in patients after acute ischemic stroke. *Oxid. Med. Cellular Longev.* doi:10.1155/2016/9761697.
29. Yildirim, E. et al. (2016). The impact of protein oxidation on sustained and white coat hypertension. *Anatol. J. Cardiol.* doi:10.14744/anatoljcardiol.2016.7174.
30. Kinoda, J. et al. (2016). Cytotoxicity of silver nanoparticle and chitin-nanofiber sheet composites caused by oxidative stress. *Nanomaterials* **6**:189.
31. Gutzkow, K. B. et al. (2016). Enhanced susceptibility of obese mice to glycidamide-induced sperm chromatin damage without increased oxidative stress. *Andrology* doi:10.1111/andr.12233.
32. Son, M. J. et al. (2016). Evaluation of the anti-fatigue effects of a traditional herbal drug, Gongjindan, under insufficient sleep conditions: study protocol for a randomised controlled trial. *Trials.* **17**:418.

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

## **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)

©2006-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.