

---

Product Manual

# OxiSelect™ AOPP Assay Kit

Catalog Number

STA-318

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 defined as an increase in the production of reactive oxygen species (ROS) due to an imbalance between antioxidant and oxidants. Advanced Oxidation Protein Products (AOPP) are uremic toxins created during oxidative stress through the reaction of chlorinated oxidants, such as chloramines and hypochlorous acid, with plasma proteins. AOPPs are structurally similar to Advanced Glycation End-Product (AGE) proteins and exert similar biological activities. AOPPs are elevated in patients with renal complications, atherosclerosis, diabetes mellitus, systemic sclerosis, as well as HIV-positive patients. Human Serum Albumin (HSA) treated with HOCl and AOPP generated *in vivo* can ignite oxidative reactions in both neutrophils and monocytes, which indicates both can be used as true mediators of inflammation. Although the mechanisms of AOPP degradation and elimination from the blood remain to be fully elucidated, it appears that the liver and spleen are mostly responsible for their isolation and removal.

The AOPP Assay has provided relevant information concerning free radical activity in many uremic associated disease states and the measurement of anti-oxidant characteristics of many compounds. The rapid and easy protocol has been modified by researchers in the evaluation of plasma and tissue samples. The AOPP-HSA concentration was determined from a Chloramine equivalence standard.

The OxiSelect™ AOPP Assay Kit offers a simple, reproducible, and consistent system for the detection of advanced oxidation protein products in plasma, lysates, and tissue homogenates. This kit includes a Chloramine standard and an AOPP Human Serum Albumin conjugate for use as a positive control. Each kit provides sufficient reagents to perform 200 tests including standard curve and unknown samples.

## **Assay Principle**

The Advanced Oxidation Protein Products (AOPP) Assay Kit is a bioassay tool for the direct quantitative measurement of AOPPs in biological samples. The unknown AOPP-containing samples or Chloramine standards are first mixed with an assay reaction initiator that begins a color development process. After a brief incubation, a stop solution is added and the samples and standards can be read with a standard colorimetric plate reader. The AOPP content in unknown samples is determined by comparison with the predetermined Chloramine standard curve.

## **Related Products**

1. STA-303: OxiSelect™ Nitrotyrosine Protein Immunoblot Kit
2. STA-305: OxiSelect™ Nitrotyrosine ELISA Kit
3. STA-308: OxiSelect™ Protein Carbonyl Immunoblot Kit
4. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit
5. STA-319: AOPP-Human Serum Albumin (AOPP-HSA)
6. STA-816: OxiSelect™ N-epsilon-(Carboxymethyl) Lysine (CML) Competitive ELISA Kit
7. STA-817: OxiSelect™ Advanced Glycation End Products (AGE) Competitive ELISA Kit

## **Kit Components**

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

1. Chloramine Standard (Part No. 231801): One 20  $\mu$ L tube of 100 mM Chloramine.
2. Chloramine Reaction Initiator (Part No. 231802): One 1.0 g bottle of powder.
3. Stop Solution (Part No. 231803): One 5 mL bottle.
4. 10X Assay Diluent (Part No. 231804): One 20 mL bottle.

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

1. AOPP-HSA Positive Control (Part No. 231805): One 100  $\mu$ L tube of 7.5 mg/mL AOPP-Human Serum Albumin at 0.14  $\mu$ mol AOPP/mg proteins.

## **Materials Not Supplied**

1. Protein samples such as purified protein, plasma, serum, cell lysates
2. Microcentrifuge and conical tubes
3. 96-well Microtiter Plate
4. Centrifuge
5. Container for preparing diluted solutions
6. Adjustable single channel micropipettes with disposable tips
7. Adjustable multichannel micropipette with disposable tips
8. Spectrophotometric microplate reader capable of reading at 340nm
9. 1X PBS

## **Storage**

Upon receipt, store the AOPP-HSA positive control at  $-20^{\circ}\text{C}$ . Store all other components at  $4^{\circ}\text{C}$ .

## **Preparation of Reagents**

- 1X Assay Diluent: Dilute the 10X Assay Diluent 1:10 with distilled or deionized water.
- Chloramine Reaction Initiator: Weigh out enough AOPP Reaction Initiator for a 200 mg/mL solution. Dissolve the powder in distilled or deionized water. Prepare only enough for the desired number of tests (eg. 100 mg dissolved in a final volume of 0.5 mL is enough to run 50 tests). It is recommended that the AOPP-HSA Positive Control be performed in duplicate each time the assay is used.  
*Note: The Chloramine Reaction Initiator solution is stable for 24-48 hours. Do not store or reuse diluted solutions.*
- AOPP-HSA Positive Control: Immediately before use, dilute an appropriate amount of the AOPP-HSA Positive Control 1:20 with 1X Assay Diluent.

## **Preparation of Standard Curve**

Dilute the Chloramine Standard 1:1000 in 1X Assay Diluent for a 100  $\mu\text{M}$  solution. Prepare a dilution series of Chloramine Standard in the concentration range of 100  $\mu\text{M}$  – 0  $\mu\text{M}$  by diluting the 100  $\mu\text{M}$  Chloramine solution in 1X Assay Diluent (Table 1). It is recommended that standards be performed in duplicate.

<b>Standard Tubes</b>	<b>Chloramine Standard (<math>\mu\text{L}</math>)</b>	<b>1X Assay Diluent (<math>\mu\text{L}</math>)</b>	<b>Chloramine Standard (<math>\mu\text{M}</math>)</b>
1	500 $\mu\text{L}$	0 $\mu\text{L}$	100
2	400 $\mu\text{L}$	100 $\mu\text{L}$	80
3	300 $\mu\text{L}$	200 $\mu\text{L}$	60
4	200 $\mu\text{L}$	300 $\mu\text{L}$	40
5	100 $\mu\text{L}$	400 $\mu\text{L}$	20
6	50 $\mu\text{L}$	450 $\mu\text{L}$	10
7	25 $\mu\text{L}$	475 $\mu\text{L}$	5
8	0 $\mu\text{L}$	500 $\mu\text{L}$	0

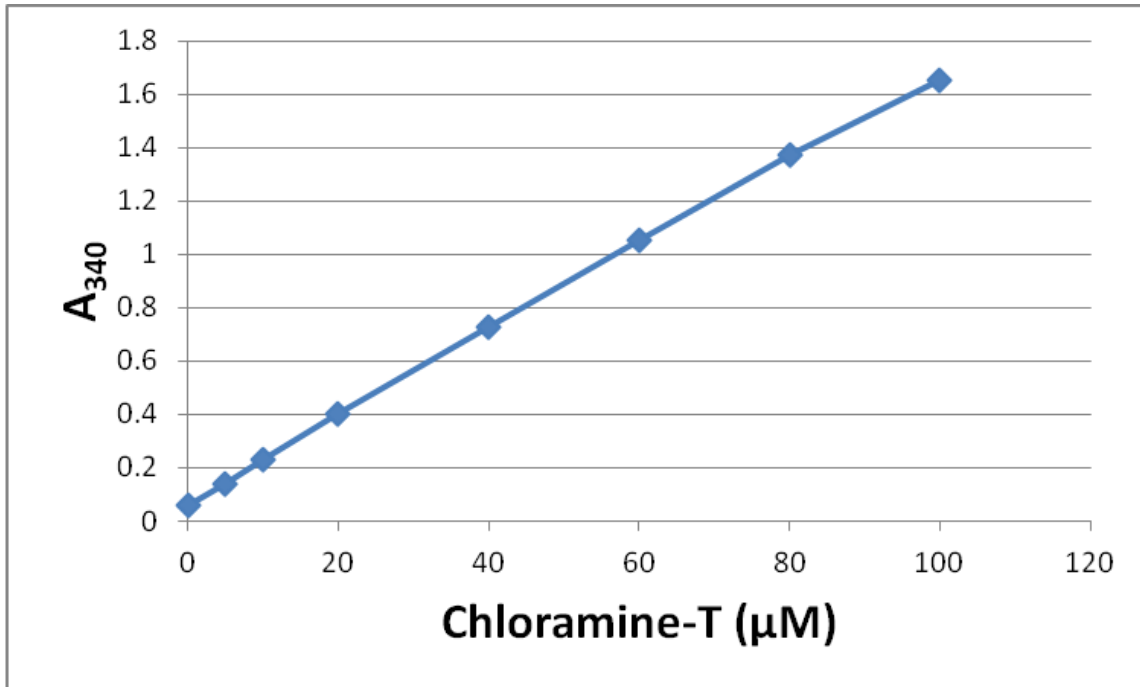
**Table 1. Preparation of Chloramine Standards**

## **Assay Protocol**

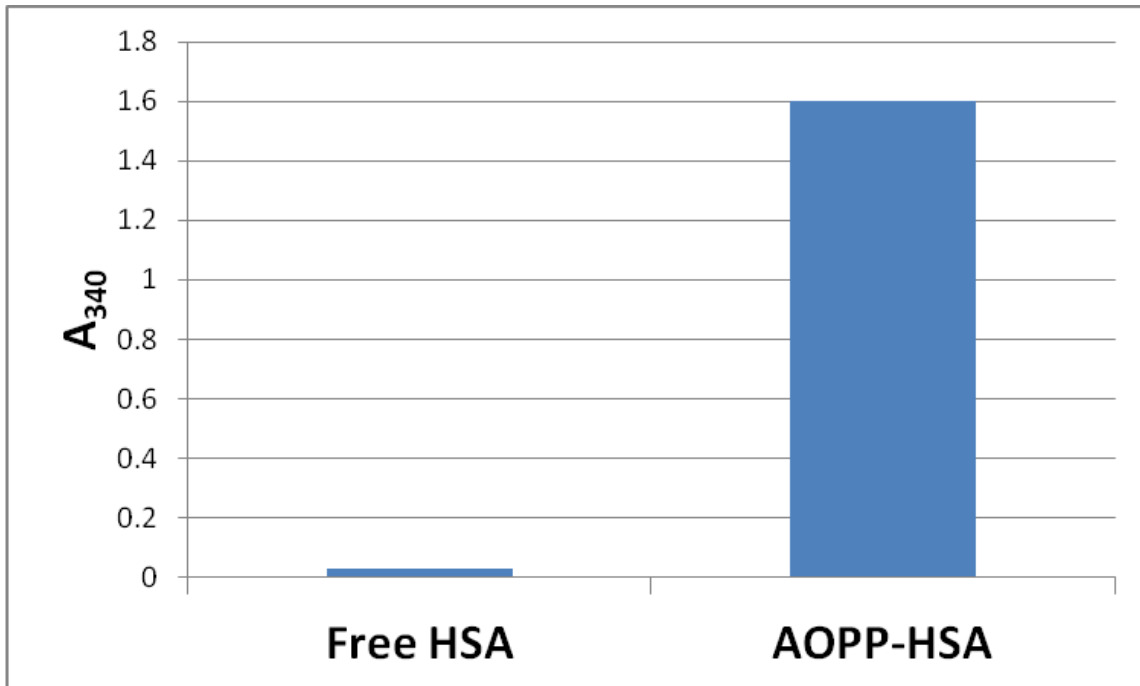
1. Prepare samples as desired. Samples such as plasma can be diluted in 1X Assay Diluent or PBS.
2. Prepare and mix all reagents thoroughly before use. Each AOPP-containing sample, the AOPP-HSA Positive Control, and Chloramine standards should be assayed in duplicate. High content AOPP samples can be further diluted for optimal analysis.
3. Add 200  $\mu\text{L}$  of samples or standards to separate wells of the microtiter plate.
4. Add 10  $\mu\text{L}$  of Chloramine Reaction Initiator to each well. Mix thoroughly and incubate on a table top rotator or shaker for 5 minutes.
5. Add 20  $\mu\text{L}$  of Stop Solution to each well. Mix thoroughly.
6. Read the absorbance of each well immediately on a spectrophotometric plate reader using 340 nm as the primary wave length. Use the 0 $\mu\text{M}$  Chloramine standard as an absorbance blank.

## **Example of Results**

The following figures demonstrate typical AOPP Assay results for the Chloramine Standard curve and the AOPP-HSA Positive Control. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1. Chloramine Standard Curve for the AOPP Assay.** The Chloramine standard curve was created as described in the Assay Protocol.



**Figure 2.** The AOPP-HSA Positive Control and untreated HSA were both prepared at a concentration of 100 µM and tested with STA-318 according to the assay protocol.

## References

1. Deschamps-Latscha B. et al. (2005) *Am. J. Kidney Dis.* 45(1): 39-47.
2. Servettaz, A. et al. (2007) *Ann. Rheum. Dis.* 66: 1202-1209.
3. Witko-Sarsat, V. et al. (1996) *Kidney International* 49: 1304-1313.

## Recent Product Citations

1. Nukala, S.B. et al. (2021). Protein network analyses of pulmonary endothelial cells in chronic thromboembolic pulmonary hypertension. *Sci Rep.* **11**(1):5583. doi: 10.1038/s41598-021-85004-z.
2. AlMarabeh, S. et al. (2020). Chronic intermittent hypoxia impairs diuretic and natriuretic responses to volume expansion in rats with preserved low-pressure baroreflex control of the kidney. *Am J Physiol Renal Physiol.* doi: 10.1152/ajprenal.00377.2020.
3. Xiao, L.L. et al. (2020). Using advanced oxidation protein products and ischaemia-modified albumin to monitor oxidative stress levels in patients with drug-induced liver injury. *Sci Rep.* **10**(1):18128. doi: 10.1038/s41598-020-75141-2.
4. Bloomer, R.J. et al. (2020). Meta- bolic Health Outcomes Following Nine Months of Mild Caloric Restriction in Male Rats Adhering To a Western or Vegan Diet. *J Altern Complement Integr Med.* **6**:097. doi: 10.24966/ACIM-7562/100097.
5. Bigazzi, R. et al. (2019). Hypertension in High School Students: Genetic and Environmental Factors: The HYGEF Study. *Hypertension.* **75**(1):71-78. doi: 10.1161/HYPERTENSIONAHA.119.13818.
6. McAllister, M.J. et al. (2019). Acute coffee ingestion with and without medium chain triglycerides decreases blood oxidative stress markers and increases ketone levels. *Can J Physiol Pharmacol.* doi: 10.1139/cjpp-2019-0458.
7. Ungurianu, A. et al. (2019). Preclinical and clinical results regarding the effects of a plant-based antidiabetic formulation versus well established antidiabetic molecules. *Pharmacol Res.* **150**:104522. doi: 10.1016/j.phrs.2019.104522.
8. Smith, C. et al. (2019). Chronic testosterone deprivation sensitizes the middle-aged rat brain to damaging effects of testosterone replacement. *Neuroendocrinology.* doi: 10.1159/000504445.
9. Gryszczyńska, B. et al. (2019). Advanced Oxidation Protein Products and Carbonylated Proteins Levels in Endovascular and Open Repair of an Abdominal Aortic Aneurysm: The Effect of Pre-, Intra-, and Postoperative Treatment. *BioMed Research International.* **2019**(7976043)1-9 pages. doi: 10.1155/2019/7976043.
10. Morsy, M.D. et al. (2019). Protective effect of combined melatonin and  $\alpha$ -tocopherol administration in spinal cord ischemia-reperfusion injury in rat. *Int. J. Morphol.* **37**(2):428-437. doi: 10.4067/S0717-95022019000200428.
11. Albatayneh, E.M. et al. (2019). Serum Oxidative-Antioxidative Status in Patients With Alkaptonuria. *J Clin Med Res.* **11**(5):337-344. doi: 10.14740/jocmr3801.
12. Shell, B. et al. (2019). Angiotensin Type 1a Receptors in the Median Preoptic Nucleus Support Intermittent Hypoxia-Induced Hypertension. *Am J Physiol Regul Integr Comp Physiol.* doi: 10.1152/ajpregu.00393.2018.
13. Bloomer, R. et al. (2018) Chronic Marijuana Smoking Does Not Negatively Impact Select Blood Oxidative Stress Biomarkers in Young, Physically Active Men and Women. *Health.* **10**:960-970. doi: 10.4236/health.2018.107071.

14. Sighinolfi, G. et al. (2018). AB0760 Advanced oxidation protein products in serum of patients with systemic sclerosis: a possible indicator of clinical evolution. *Annals of the Rheumatic Diseases*. **77**:1516. doi: 10.1136/annrheumdis-2018-eular.6809.
15. Wilson, E.N. et al. (2018). Chronic intermittent hypoxia induces hormonal and male sexual behavioral changes: Hypoxia as an advancer of aging. *Physiol Behav*. **189**:64-73. doi: 10.1016/j.physbeh.2018.03.007.
16. Gradinaru, D. et al. (2018). Insulin-Leptin Axis, Cardiometabolic Risk and Oxidative Stress in Elderly with Metabolic Syndrome. *Exp Clin Endocrinol Diabetes*. doi: 10.1055/s-0043-123825.
17. Hányšová, S. et al. (2017). Elevated plasma levels of advanced oxidation protein products in Slovak multiple sclerosis patients: possible association with different disability states. *Acta Nerv Super Rediviva*. **59**(2): 45–50.
18. Sun, S. et al. (2018). Advanced oxidation protein products induce S-phase arrest of hepatocytes via the ROS-dependent,  $\beta$ -catenin-CDK2-mediated pathway. *Redox Biol*. **14**:338-353. doi: 10.1016/j.redox.2017.09.011.
19. Snyder, B. et al. (2017). Chronic intermittent hypoxia induces oxidative stress and inflammation in brain regions associated with early-stage neurodegeneration. *Physiol. Rep*. doi:10.14814/phy2.13258.
20. Budzyń, M. et al. (2017). The Association of Serum Thrombomodulin with Endothelial Injuring Factors in Abdominal Aortic Aneurysm. *Hindawi BioMed Research International*. doi: 10.1155/2017/2791082.
21. Gradinaru, D. et al. (2016). Adiponectin: possible link between metabolic stress and oxidative stress in the elderly. *Aging Clin Exp Res*. doi:10.1007/s40520-016-0629-z.
22. Wan, X. et al. (2016). SIRT1-PGC1 $\alpha$ -NF $\kappa$ B pathway of oxidative and inflammatory stress during *Trypanosoma cruzi* infection: benefits of SIRT1-targeted therapy in improving heart function in Chagas disease. *PLoS Pathog*. doi: 10.1371/journal.ppat.1005954.
23. Crone, L. B. et al. (2016). Impact of meal ingestion rate and caffeine coingestion on postprandial lipemia and oxidative stress following high-fat meal consumption. *J Caffeine Res*. doi:10.1089/jcr.2016.0004.
24. Huemer, M. et al. (2016). Clinical phenotype, biochemical profile, and treatment in 19 patients with arginase 1 deficiency. *J Inherit Metab Dis*. doi:10.1007/s10545-016-9928-y.
25. Martins, L. S. et al. (2015). Advanced Glycation Endproducts (AGE) evolution after pancreas-kidney transplantation: plasmatic and cutaneous assessments. *Oxid Med Cell Longev*. 2189582.
26. Šarić, B. et al. (2015). Oxidative stress impact on growth hormone secretion in the eye. *Croat Med J*. **56**:326-333.
27. Bloomer, R. J. et al. (2015). Comparison of a restricted and unrestricted vegan diet plan with a restricted omnivorous diet plan on health-specific measures. *Healthcare*. **3**:544-555.
28. Jung, E. et al. (2015). Gemigliptin improves renal function and attenuates podocyte injury in mice with diabetic nephropathy. *Eur J Pharmacol*. doi: 10.1016/j.ejphar.2015.04.055.
29. Thurmond, P. et al. (2015). Structural modifications of the prostate in hypoxia, oxidative stress, and chronic ischemia. *Korean J Urol*. **56**:187-196.
30. Xie, Z. X. et al. (2014). Effect of GABA on oxidative stress in the skeletal muscles and plasma free amino acids in mice fed high-fat diet. *J Anim Physiol Anim Nutr (Berl)*. doi: 10.1111/jpn.12254.

## **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)

©2007-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.