

---

Product Manual

# OxiSelect™ Ferric Reducing Antioxidant Power (FRAP) Assay Kit

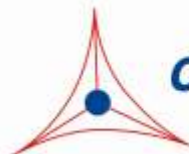
Catalog Number

STA-859

200 assays

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

---



**CELL BIOLABS, INC.**

*Creating Solutions for Life Science Research*

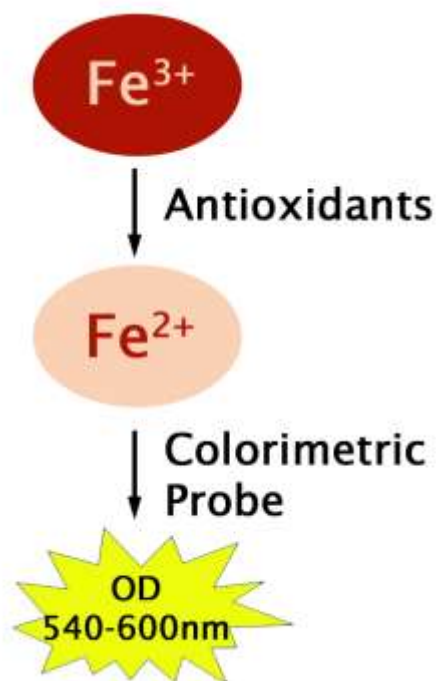
## **Introduction**

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 inflammation and 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 are molecules that contain one or more free electrons that can be donated to stabilize ROS. Antioxidants 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 neutralized and eliminated from the cells by this variety of antioxidant defense mechanisms.

Cell Biolabs' OxiSelect™ Ferric Reducing Antioxidant Power (FRAP) Assay Kit is a quantitative assay for measuring the antioxidant potential within various samples such as serum, plasma, urine, saliva, tears, tissue homogenates, cell extracts, and purified food or drug extracts. The assay is based on the highly-cited work of Benzie and Strain (ref. 1 and 2) in which iron reacts with a colorimetric probe to produce a blue product. The reaction is driven by the electron donating reducing power of antioxidants. The kit employs a ferrous iron standard which allows the user to determine the antioxidant content present within their sample. Each kit provides sufficient reagents to perform up to 200 assays, including standard curve and unknown samples. The FRAP assay offers a simple and efficient analytical method for assessing age, disease, diet, or other physiological changes to antioxidant status.

## **Assay Principle**

The OxiSelect™ Ferric Reducing Antioxidant Power (FRAP) Assay Kit is a quantitative assay for measuring the antioxidant potential within a sample. Ferric iron ( $\text{Fe}^{3+}$ ) is initially reduced by electron-donating antioxidants present within the sample to its ferrous form ( $\text{Fe}^{2+}$ ). The iron-colorimetric probe complex develops a dark blue color product upon reduction which can be measured at 540-600 nm (see Figure 1). Samples can be compared to the iron standard for determining antioxidant potential. This assay is analytically sensitive to approximately 4  $\mu\text{M}$  of  $\text{Fe}^{2+}$  iron equivalents, or a FRAP value of 2.



**Figure 1. Assay Principle.**

### **Related Products**

1. STA-312: OxiSelect™ Total Glutathione Assay Kit
2. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
3. STA-340: OxiSelect™ Superoxide Dismutase Activity Assay
4. STA-341: OxiSelect™ Catalase Activity Assay Kit
5. STA-342: OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence)
6. STA-345: OxiSelect™ ORAC Activity Assay
7. STA-347: OxiSelect™ *In Vitro* ROS/RNS Assay Kit (Green Fluorescence)
8. STA-360: OxiSelect™ Total Antioxidant Capacity (TAC) Assay Kit
9. STA-802: OxiSelect™ *In Vitro* Nitric Oxide (Nitrite / Nitrate) Assay kit
10. STA-803: OxiSelect™ Myeloperoxidase Activity Assay Kit (Colorimetric)
11. STA-812: OxiSelect™ Glutathione Reductase Assay Kit
12. STA-860: OxiSelect™ Ascorbic Acid Assay Kit (FRASC) (Colorimetric)

### **Kit Components**

1. Colorimetric Probe (Part No. 286001): Two 1 mL tubes
2. Iron Chloride Solution (Part No. 286002): Two 1 mL tubes
3. Iron (II) Standard (Part No. 285901): One 500 mg amber tube of crystals
4. Assay Buffer (5X) (Part No. 286004): Two 25 mL bottles

## **Materials Not Supplied**

1. 96-well microtiter plates
2. 10 kDa MWCO centrifugal filter (for high protein content samples)
3. Ethanol or other organic solvent for lipid-based samples
4. 10  $\mu$ L to 1000  $\mu$ L adjustable single channel micropipettes with disposable tips
5. 50  $\mu$ L to 300  $\mu$ L adjustable multichannel micropipette with disposable tips
6. Conical tubes, microcentrifuge tubes, and bottles for sample and buffer preparation
7. Centrifuge and/or microfuge
8. Sonicator or tissue homogenizer
9. Multichannel micropipette reservoirs
10. Microplate reader capable of reading absorbance between 540-600 nm

## **Storage**

Upon receipt, store the kit components at 4°C.

## **Preparation of Reagents**

- 1X Assay Buffer: Prepare 1X Assay Buffer by adding deionized water to the 5X Assay Buffer (e.g. add 3 mL of 5X Assay Buffer to 12 mL of deionized water). Mix thoroughly until homogeneous. Use this buffer for preparing kit reagents and within the assay. Store at 4°C when not in use.
- Reaction Reagent: Prepare the Reaction Reagent just before use and prepare only enough for immediate applications. Dilute the Colorimetric Probe 1:10 and the Iron Chloride Solution 1:10 in 1X Assay Buffer (e.g. For 50 assays, add 500  $\mu$ L Colorimetric Probe and 500  $\mu$ L Iron Chloride Solution to 4 mL 1X Assay Buffer for 5 mL total). Vortex thoroughly.

## **Preparation of Samples**

These preparation protocols are intended as a guide for preparing unknown samples. The user may need to adjust the sample treatment accordingly. All samples should be assayed immediately or stored for up to 2 months at -80°C. A trial assay with a representative test sample should be performed to determine the sample compatibility with the dynamic range of the standard curve. High levels of interfering substances may cause variations in results. Samples may be diluted in 1X Assay Buffer as necessary before testing. Run proper controls as necessary. Always run a standard curve with samples.

- Serum\*: Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C.
- Plasma\*: Collect blood sample and add to a blood collection tube containing heparin as the anticoagulant. Centrifuge at 3,000 rpm for 10-15 minutes at 4°C. Remove the upper yellow plasma supernatant layer without disturbing the white buffy coat (leukocytes). Samples should be tested immediately or frozen at -80°C.

*\*Note: Hemolyzed plasma or serum should be avoided. Heparinized plasma is recommended over EDTA plasma.*

- Cell lysates: Lyse 1-2 x 10<sup>6</sup> cells/mL by sonication or multiple freeze-thaw cycles in 4 volumes of cold 1X Assay Buffer. Centrifuge at 12,000 rpm for 15 minutes at 4°C and remove insoluble cell material. A high concentration of protein may interfere with the assay. In this case, filter the sample with a 10kDa MWCO centrifugal filter before assaying (to reduce protein interference and turbidity). Test samples immediately or store at -80°C.
- Tissue lysates/homogenates: Homogenize/sonicate approximately 10 mg of tissue in 1-2 mL cold 1X Assay Buffer. Centrifuge the homogenate at 12,000 rpm for 15 minutes at 4°C and collect the supernatant. A high concentration of protein may interfere with the assay. In this case, filter the sample with a 10kDa MWCO centrifugal filter before assaying (to reduce protein interference and turbidity). Test samples immediately or store at -80°C.
- Food extracts: Wash uncooked food (e.g. fruits or vegetables) in water and homogenize 5-10g (fresh wet weight) in 100 mL distilled water for 30 seconds. Filter homogenate and prepare dilutions in deionized water, 70% acetone, or 50% ethanol as necessary.

### **Preparation of Standard Curve**

*Note: The Iron (II) Standards should be prepared immediately prior to running the assay.*

Prepare Iron (II) Standards fresh by weighing out the iron standard crystals for a 10 mg/mL solution in deionized water. This 10 mg/mL solution is equivalent to a concentration of 36 mM. Next, dilute this 36 mM stock to 1 mM in deionized water (e.g. add 125 µL of the 36 mM iron(II) solution to 4.375 mL deionized water). This is equivalent to a 1 mM or 1000 µM solution. Use this solution to prepare a series of standards according to Table 1 below. Prepare standards immediately prior to each assay performed. Vortex tubes thoroughly. Do not store or reuse the standard preparations.

<b>Standard Tubes</b>	<b>Iron (II) 1000 µM (µL)</b>	<b>Deionized Water (µL)</b>	<b>Iron (II) Concentration (µM)</b>
1	500	500	500
2	500 of tube #1	500	250
3	500 of tube #2	500	125
4	500 of tube #3	500	62.5
5	500 of tube #4	500	31.3
6	500 of tube #5	500	15.6
7	500 of tube #6	500	7.8
8	500 of tube #7	500	3.9
9	0	500	0

**Table 1. Preparation of Iron(II) Standards.**

*Note: The Iron (II) Standard solutions are oxidized over time, yielding a yellow precipitate. Discard all iron solutions after use. Do not use iron solutions that contain precipitate.*

## Assay Protocol

1. Prepare and mix all reagents thoroughly before use. Each standard, sample and control should be assayed in duplicate or triplicate.
2. Add 100  $\mu\text{L}$  of each standard, unknown sample or control to a 96-well plate.
3. Add 100  $\mu\text{L}$  of the Reaction Reagent to all wells and mix by pipetting or with a horizontal shaker. Immediately time the reaction after adding the reagent and incubate 10 minutes at room temperature.

*Note: Reaction times may vary for some antioxidants. The assay reaction time can be extended, but a longer reaction time will result in higher background.*

4. Immediately read absorbance of each well on a microplate reader using 540-600 nm as the primary wave length.

## Calculation of Results

1. Determine the average absorbance values for each sample, control, and standard.
2. Calculate the net OD by subtracting the zero standard value from samples and standards. This is the background correction.
3. Graph the standard curve (see Figure 2).
4. Compare the net OD of each sample to the standard curve to determine the quantity of antioxidant potential, as  $\mu\text{M Fe}^{2+}$  iron equivalents (FRAP value), present in the sample. Only use values within the range of the standard curve.

## Example of Results

The following figures demonstrate typical FRAP Assay results at 540 nm. One should use the data below for reference only. This data should not be used to interpret actual results.

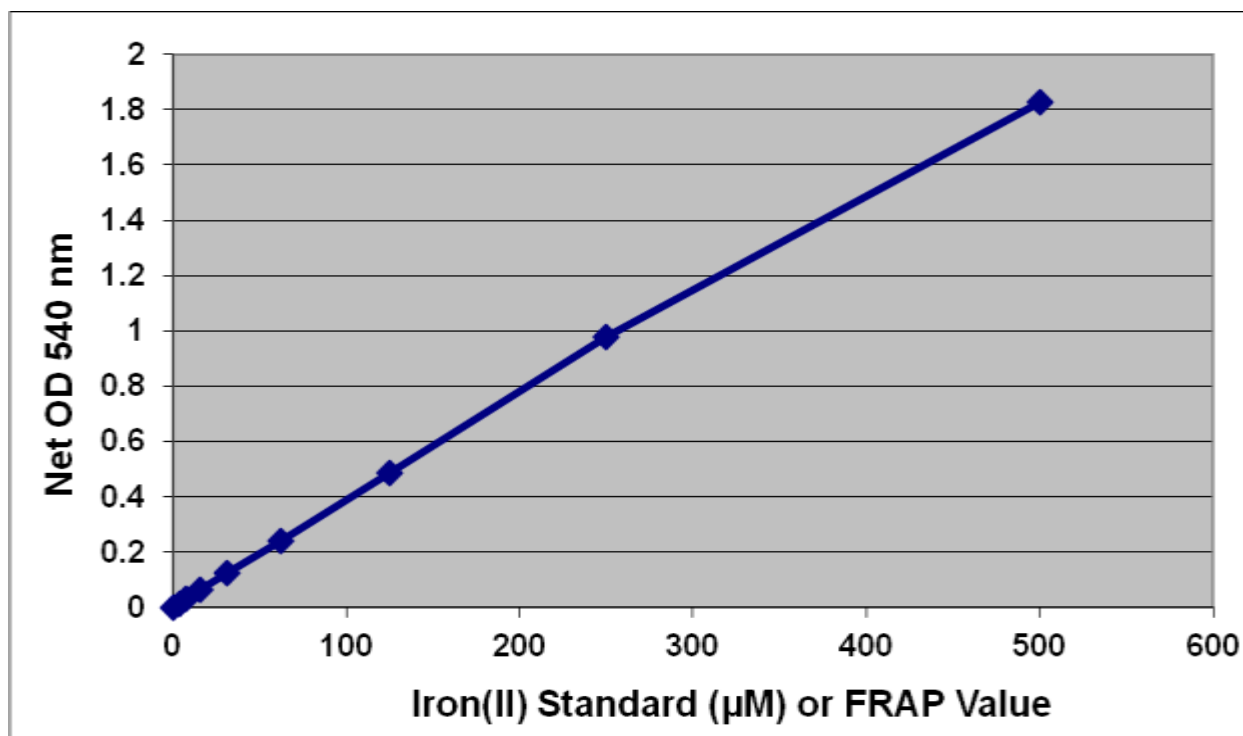
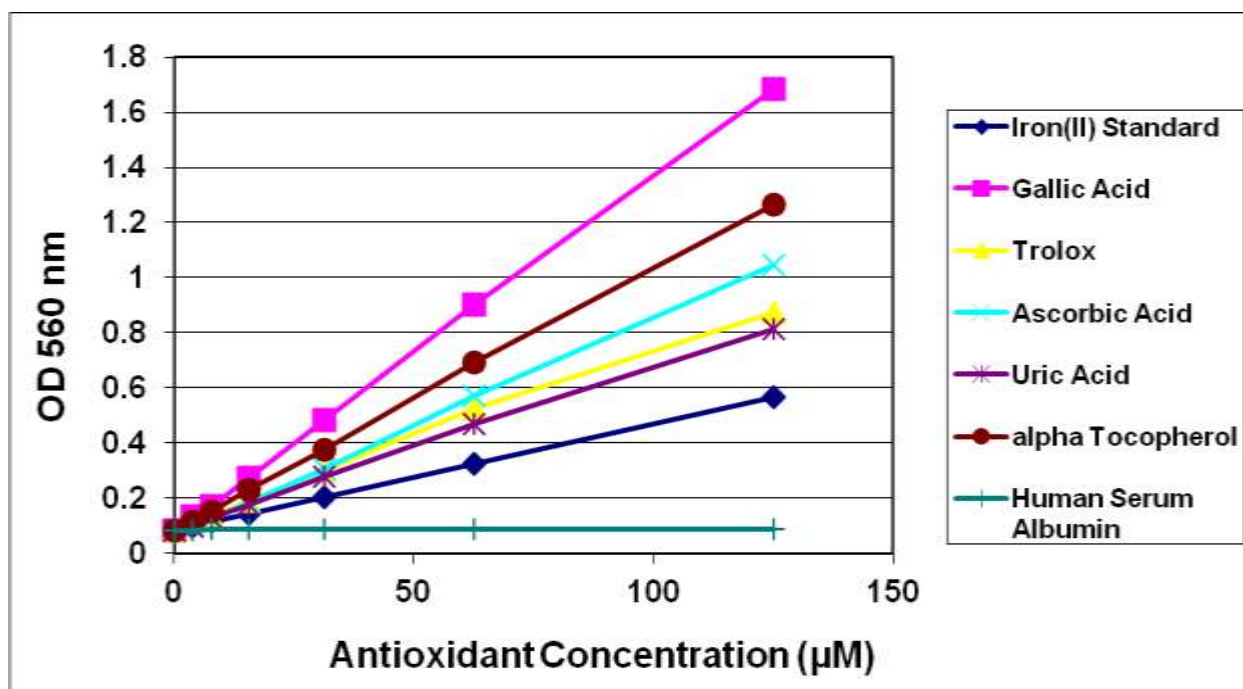


Figure 2. Iron (II) Standard Curve.



**Figure 3. FRAP Assay results for various antioxidants.**

### References

1. Benzie, I.F.F., et. al. *Redox Report* (1997) 3: 233-238.
2. Benzie, I.F.F., et. al. *Methods Enzymol.* (1999) 299: 15-23.
3. Jenab, M., et. al. *Cancer Epidemiol. Bio. Prev.* (2005) 14(7): 1837-1840.
4. Karlsen, A., et al. *Eur. J. Clin. Nut.* (2007) 61: 1233-1236.
5. Siu, P. M., et. al. *Exp Physiol.* (2011) 96.9: 889-906.

### Recent Product Citations

1. DeJulius, C.R. et al. (2021). Optimizing an Antioxidant TEMPO Copolymer for Reactive Oxygen Species Scavenging and Anti-Inflammatory Effects in Vivo. *Bioconjug Chem.* doi: 10.1021/acs.bioconjugchem.1c00081.
2. He, H. et al. (2021). Vinegar/Tetramethylpyrazine Induces Nutritional Preconditioning Protecting the Myocardium Mediated by VDAC1. *Oxid Med Cell Longev.* doi: 10.1155/2021/6670088.
3. Yeh, W.J. et al. (2020). *Hylocereus polyrhizus* Peel Extract Retards Alcoholic Liver Disease Progression by Modulating Oxidative Stress and Inflammatory Responses in C57BL/6 Mice. *Nutrients.* **12**(12):3884. doi: 10.3390/nu12123884.
4. Nakov, G. et al. (2020). Effect of grape pomace powder addition on chemical, nutritional and technological properties of cakes. *LWT.* doi: 10.1016/j.lwt.2020.109950.
5. Belmans, N. et al. (2020). Quantification of DNA Double Strand Breaks and Oxidation Response in Children and Adults Undergoing Dental CBCT Scan. *Sci Rep.* **10**(1):2113. doi: 10.1038/s41598-020-58746-5.
6. Yang, B. et al. (2019). Tetramethylpyrazine Attenuates the Endotheliotoxicity and the Mitochondrial Dysfunction by Doxorubicin via 14-3-3γ/Bcl-2. *Oxid Med Cell Longev.* doi: 10.1155/2019/5820415.

7. Balgoon, M.J. et al. (2019). Assessment of the Protective Effect of *Lepidium sativum* against Aluminum-Induced Liver and Kidney Effects in Albino Rat. *BioMed Research International*. doi: 10.1155/2019/4516730.
8. Luo, Y. et al. (2019). Nutritional preconditioning induced by astragaloside IV on isolated hearts and cardiomyocytes against myocardial ischemia injury via improving Bcl-2-mediated mitochondrial function. *Chem Biol Interact*. pii: S0009-2797(19)30320-5. doi: 10.1016/j.cbi.2019.06.036.
9. Belmans, N. et al. (2019). Method validation to assess in vivo cellular and subcellular changes in buccal mucosa cells and saliva following CBCT examinations. *Dentomaxillofac Radiol*. 20180428. doi: 10.1259/dmfr.20180428.
10. Innes, S.N. et al. (2019). Elevated air humidity increases UV mediated leaf and DNA damage in pea (*Pisum sativum*) due to reduced flavonoid content and antioxidant power. *Photochem Photobiol Sci*. **18**(2):387-399. doi: 10.1039/c8pp00401c.
11. Hosokawa, Y. et al. (2016). Ascorbic acid does not reduce the anticancer effect of radiotherapy. *Biomed. Rep*. doi:10.3892/br.2016.819.

## **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)

©2015-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.