Product Manual

OxiSelect[™] 8-iso-Prostaglandin F2α ELISA Kit

Catalog Numbers

STA-337	96 assays
STA-337-5	5 x 96 assays

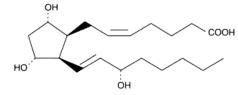
FOR RESEARCH USE ONLY Not for use in diagnostic procedures



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 isoprostanes. The isoprostanes are a type of eicosanoids produced non-enzymatically through the oxygen radical induced peroxidation of tissue phospholipids and lipoproteins. Isoprostanes are prostaglandin-like compounds that appear in normal plasma and urine samples, but are elevated by oxidative stress in tissue, plasma, and urine.

8-iso-Prostaglandin F2 α (also known as 8-epi-PGF2 α , 8-isoprostane, or 15-isoprostane F2t), is an isoprostane that has been shown to be useful for the assessment of oxidative stress *in vivo*. It is produced in membrane phospholipids from non-cyclooxygenase and cyclooxygenase peroxidation pathways derived from arachidonic acid. 8-iso-Prostaglandin F2 α (8-iso-PGF2 α) is a potent vasoconstrictor, a mutagen in 3T3 cells as well as vascular smooth muscle cells, and also a possible pathophysiological mediator that can alter membrane integrity. It has been implicated in atherogenesis and elevated levels are associated with hepatorenal syndrome, rheumatoid arthritis, carcinogenesis, as well as atherosclerosis. 8-iso-PGF2 α circulates in the plasma and is excreted in the urine. 8-iso PGF2 α circulates as an esterified LDL Phospholipid and as a free acid. Normal human plasma and urine 8-iso PGF2 α is about 40-100 pg/mL and about 190 pg/mg of creatinine respectively. Methods for determining total 8-iso PGF2 α usually require alkaline hydrolysis of 8-iso PGF2 α esters from tissues followed by extractions, phase separations and thin layer chromatography.



8-iso-Prostaglandin F2a (8-iso-PGF2a)

The OxiSelectTM 8-iso-Prostaglandin F2 α ELISA Kit is an enzyme immunoassay developed for rapid detection and quantification of 8-iso-PGF2 α . The quantity of 8-iso-PGF2 α in samples is determined by comparing its absorbance with that of a known 8-iso-PGF2 α standard curve. Each kit provides sufficient reagents to perform up to 96 assays, including the standard curve and unknown samples.

Assay Principle

Cell Biolabs' 8-iso-PGF2 α kit is a competitive enzyme-linked immunoassay (ELISA) for determining levels of 8-iso-PGF2 α in a variety of biological samples such as plasma, urine, serum, or tissue extracts. An antibody to 8-iso-PGF2 α is incubated in pre-coated microtiter plate wells. Upon washing, 8-iso-PGF2 α standards or treated samples are mixed with an 8-iso-PGF2 α -HRP conjugate and added simultaneously to the wells. The unconjugated, or free 8-iso-PGF2 α and 8-iso-PGF2 α -HRP conjugate compete for binding to the antibody bound to the plate. After this brief incubation and wash, a substrate to the HRP is added. The HRP activity results in color development that is directly proportional to the amount of 8-iso-PGF2 α in the samples or standards. The 8-iso-PGF2 α content in an unknown sample is determined by comparing with the known predetermined standard curve. Please read the complete kit insert prior to performing the assay.



Related Products

- 1. STA-320: OxiSelectTM Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
- 2. STA-325: OxiSelect[™] Oxidative RNA Damage ELISA Kit (8-OHG Quantitation)
- 3. STA-330: OxiSelect[™] TBARS Assay Kit (MDA Quantitation)
- 4. STA-344: OxiSelect[™] Hydrogen Peroxide/Peroxidase Assay Kit
- 5. STA-347: OxiSelectTM In Vitro ROS/RNS Assay Kit (Green Fluorescence)
- 6. STA-816: OxiSelectTM N-epsilon-(Carboxymethyl) Lysine (CML) Competitive ELISA Kit
- 7. STA-817: OxiSelect[™] Advanced Glycation End Products (AGE) Competitive ELISA Kit
- 8. STA-832: OxiSelect[™] MDA Competitive ELISA Kit
- 9. STA-838: OxiSelect[™] HNE Adduct Competitive ELISA Kit

Kit Components

- 1. Goat Anti-Rabbit Antibody Coated Plate (Part No. 250001): One 96-well strip plate.
- 2. <u>Anti-8-iso-PGF2α Antibody</u> (Part No. 233701): One 20 μL tube of anti-8-iso-PGF2α rabbit IgG.
- 3. <u>Sample Diluent</u> (Part No. 233702): One 50 mL bottle.
- 4. <u>Neutralization Solution</u> (Part No. 233705): One 20 mL bottle.
- 5. <u>10X Wash Buffer</u> (Part No. 310806): One 100 mL bottle.
- 6. <u>Substrate Solution</u> (Part No. 310807): One 12 mL amber bottle.
- 7. <u>Stop Solution</u> (Part. No. 310808): One 12 mL bottle.
- 8. <u>8-iso-PGF2α Standard</u> (Part No. 233703): One 25 μL tube of 200 μg/mL 8-iso-PGF2α in DMSO.
- 9. <u>8-iso-PGF2α-HRP Conjugate</u> (Part No. 233704): One 70 μL tube of 8-iso-PGF2α-HRP conjugate.

Materials Not Supplied

- 1. Protein samples such as purified protein, plasma, serum, cell lysate
- 2. Deionized water
- 3. $5 \,\mu$ L to 1000 μ L adjustable single channel precision micropipettes with disposable tips
- 4. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
- 5. Bottles, flasks, and conical or microtubes necessary for reagent preparation
- 6. Reagents and materials necessary for sample extraction and purification
- 7. Multichannel micropipette reservoir
- 8. Plate orbital shaker or rotator
- 9. Microplate reader capable of reading at 450 nm (620 nm as optional reference wave length)



Storage

Upon receipt, store the Anti-8-iso-PGF2α Antibody, 8-iso-PGF2α-HRP Conjugate, and 8-iso-PGF2α Standard at -20°C. Make aliquots as necessary to avoid freeze/thaw cycles. Store all other kit components at 4°C. Any partial or unused components should return to their proper storage temperatures.

Safety Considerations

- Some kit components contain azide, which can react with copper or lead piping. Flush with large volumes of water when disposing of reagents.
- Some kit reagents are caustic or hazardous and should be handled accordingly.

Preparation of Reagents

- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.
- Anti-8-iso-PGF2α Antibody: Immediately before use, dilute the Anti-8-iso-PGF2α Antibody 1:1000 with Sample Diluent.
- 8-iso-PGF2α-HRP Conjugate: Immediately before use, dilute the conjugate 1:80 with Sample Diluent. Only prepare enough of the diluted conjugate for the number of wells immediately used.
- Substrate Solution: Prior to use, warm the Substrate Solution to room temperature.

Note: Do not store diluted Anti-8-iso-PGF2a Antibody, 8-iso-PGF2a-HRP Conjugate, or 8-iso-PGF2a Standard solutions.

Preparation of Samples

Hydrolysis of lipoprotein or phospholipid coupled 8-iso-Prostaglandin F2 α (8-iso-PGF2 α) is required to measure both free and esterified isoprostane. To hydrolyze this ester bond, the sample is usually treated with 2N NaOH at 45°C for 2 hours.

- Serum, plasma, tissue lysate samples: Use 1 part of 10N NaOH for every 4 parts of liquid sample. After incubation at 45°C for 2 hours, add 100 μL of concentrated (10N) HCl per 500 μL of hydrolyzed sample. The sample could turn milky after this addition. Centrifuge the samples for 5 minutes at 12,000 rpm in a microcentrifuge. The clear supernatant can be used in the assay or stored at -20°C or below for future use. Before assaying, check to be sure each neutralized sample is in the pH range of 6-8. If it is not, adjust the pH to this range by adding 100 μL of the sample to 100 μL of the provided Neutralization Solution.
- Urine samples: Acid hydrolysis of urine samples is necessary to break the bonds which hold lipid and non-lipid components together prior to ELISA. Urine sample is acidified to pH 3.0 by adding 1/10 volume of 1N HCl (Example: Add 100 μ L of 1N HCl to 1 mL of urine sample). Acidified urine sample should be further diluted in PBS or Sample Diluent 1:4 to 1:8 before ELISA.



Preparation of 8-iso-PGF2α Standards

1. Prepare fresh standards by diluting the 8-iso-PGF2 α Standard from 200 μ g/mL to 0.2 μ g/mL in Sample Diluent for a 1:1000 final dilution. (Example: Add 5 μ L of 8-iso-PGF2 α Standard stock tube to 4.995 mL of Sample Diluent)

Standard Tubes	8-iso-PGF2α Standard (μL)	Sample Diluent (µL)	8-iso-PGF2α Standard (pg/mL)
1	5 µL of Standard Stock	4995 μL	200,000
2	250 μL of Tube #1	750 μL	50,000
3	250 μL of Tube #2	750 μL	12,500
4	250 μL of Tube #3	750 μL	3,125
5	250 μL of Tube #4	750 μL	781
6	250 μL of Tube #5	750 µL	195
7	250 μL of Tube #6	750 μL	49
8	0 µL	200 µL	0

2. Prepare a series of the remaining 8-iso-PGF2 α standards according to Table 1.

Table 1. Preparation of 8-iso-PGF2α Standard Curve.

Note: Do not store diluted 8-iso-PGF2a Standard solutions.

Assay Protocol

Note: Each 8-iso-PGF2 α Standard and unknown samples should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

- Add 100 μL of the diluted Anti-8-iso-PGF2α Antibody to the Goat Anti-Rabbit Antibody Coated Plate. Incubate 1 hour at 25°C on an orbital shaker.
- Remove the antibody solution from the wells. Wash wells 5 times with 300 µL 1X Wash Buffer per well. After the last wash, empty the wells and tap microwell plate on absorbent pad or paper towel to remove excess wash solution.

Note: Thorough washing is necessary to remove all of the azide present in the antibody solution.

- 3. Combine 55 μ L of the 8-iso-PGF2 α standard or sample and 55 μ L of 8-iso-PGF2 α -HRP conjugate in a microtube and mix thoroughly. Transfer 100 μ L of the combined solution per well. A well containing Sample Diluent can be used as a control. Incubate 1 hour at 25°C on an orbital shaker.
- Remove the combined solution from the wells. Wash 5 times with 300 μL of 1X Wash Buffer per well. After the last wash, empty wells and tap microwell plate on absorbent pad or paper towel to remove excess wash solution.



- 5. Add 100 μ L of Substrate Solution to each well. Incubate at room temperature for 10-30 minutes on an orbital shaker.
- 6. Stop the enzyme reaction by adding 100 μ L of Stop Solution to each well. Results should be read immediately (color will fade over time).
- 7. Read absorbance of each well on a microplate reader using 450 nm as the primary wave length.

Example of Results

The following figures demonstrate typical 8-iso-PGF2 α results. One should use the data below for reference only. This data should not be used to interpret actual results.

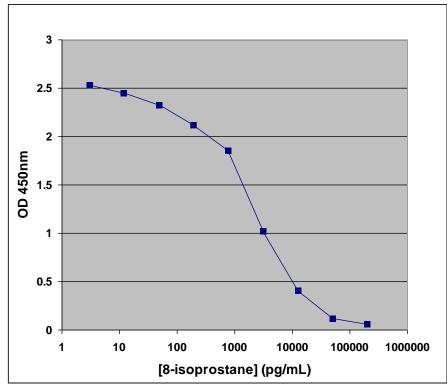


Figure 1: 8-iso-PGF2a ELISA Standard Curve.



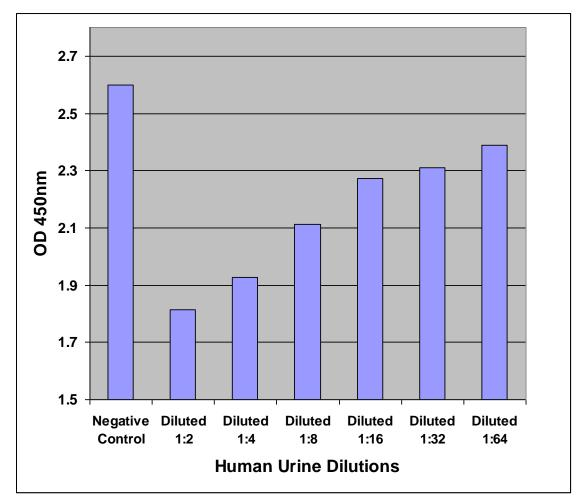


Figure 2: Dilutions of Human Urine tested with 8-iso-PGF2a ELISA.

Cross reactivity of 8-iso-Prostaglandin F2a ELISA Kit

<u>Compounds</u>	Cross Reactivity
8-iso-PGF2α	100%
PGF1a	4.6%
PGF2a	1.85%
PGE1	0.19%
TXB2	0.023%
PGB1	0.02%
PGE3	0.012%
6-keto-PGF1α	0.008%
13,14-dihydro-15-keto-PGF2α	0.008%
6,15-keto-13,14-dihydro-PGF1α	0.005%
8-iso-PGE1	<0.001%
PGA2	<0.001%
PGJ2	<0.001%



References

- 1. Banerjee, M., Kang, K.H., Morrow, J.D., et al. (1992) Am. J. Physiol. 263: H660-H663.
- 2. Morrow, J.D., Hill, K.E., Burk, R.F., et al. (1990) Proc. Natl. Acad. Sci. USA. 87: 9383-9387.
- 3. Morrow, J.D., Harris, T.M., Roberts, L.J. (1990) Anal. Biochem. 184: 1-10.
- 4. Vacchiano, C.A., and Tempel, G.E. (1994) J. Appl. Physiol. 77: 2912-2917.
- 5. Wang, Z., Ciabattoni, G., Cre'minon, C., et al. (1995) Pharmacol. Exp. Ther. 275: 94-100.

Recent Product Citations

- 1. Kopacz, A. et al. (2023). Co-administration of angiotensin II and simvastatin triggers kidney injury upon heme oxygenase-1 deficiency. *Free Radic Biol Med.* **205**:188-201. doi: 10.1016/j.freeradbiomed.2023.05.018.
- 2. Zhang, C. et al. (2023). Liver fibrosis is a common pathological change in the liver of dairy cows with fatty liver. *J Dairy Sci.* **106**(4):2700-2715. doi: 10.3168/jds.2022-22021.
- Rodriguez-Pérez, M.D. et al. (2023). The Effect of the Extra Virgin Olive Oil Minor Phenolic Compound 3',4'-Dihydroxyphenylglycol in Experimental Diabetic Kidney Disease. *Nutrients*. 15(2):377. doi: 10.3390/nu15020377.
- 4. Kumar, P. et al. (2022). Supplementing Glycine and N-Acetylcysteine (GlyNAC) in Older Adults Improves Glutathione Deficiency, Oxidative Stress, Mitochondrial Dysfunction, Inflammation, Physical Function, and Aging Hallmarks: A Randomized Clinical Trial. *J Gerontol A Biol Sci Med Sci.* doi: 10.1093/gerona/glac135.
- 5. Cao, N. et al. (2022). The Activated AMPK/mTORC2 Signaling Pathway Associated with Oxidative Stress in Seminal Plasma Contributes to Idiopathic Asthenozoospermia. *Oxid Med Cell Longev*. doi: 10.1155/2022/4240490.
- Tsunenaga, M. et al. (2022). Modulating effects of oral administration of Lycii Fructus extracts on UVB-induced skin erythema: A Randomized, placebo-controlled study. *Biomed Rep.* 17(1):62. doi: 10.3892/br.2022.1545.
- 7. Hoferichter, F. & Raufelder, D. (2022). Biophysiological stress markers relate differently to grit and school engagement among lower- and higher-track secondary school students. *Br J Educ Psychol*. doi: 10.1111/bjep.12514.
- 8. Makris, K.C. et al. (2022). Oxidative stress of glyphosate, AMPA and metabolites of pyrethroids and chlorpyrifos pesticides among primary school children in Cyprus. *Environ Res.* **212**(Pt B):113316. doi: 10.1016/j.envres.2022.113316.
- Maciejczyk, M. et al. (2022). α-Lipoic Acid Strengthens the Antioxidant Barrier and Reduces Oxidative, Nitrosative, and Glycative Damage, as well as Inhibits Inflammation and Apoptosis in the Hypothalamus but Not in the Cerebral Cortex of Insulin-Resistant Rats. Oxid Med Cell Longev. doi: 10.1155/2022/7450514.
- 10. Cañizo Vázquez, D. et al. (2022). Oxidative Stress and Indicators of Brain Damage Following Pediatric Heart Surgery. *Antioxidants (Basel)*. **11**(3):489. doi: 10.3390/antiox11030489.
- 11. Konstantinou, C. et al. (2022). Use of metabolomics in refining the effect of an organic food intervention on biomarkers of exposure to pesticides and biomarkers of oxidative damage in primary school children in Cyprus: A cluster-randomized cross-over trial. *Environ Int*. doi: 10.1016/j.envint.2021.107008.
- 12. Kumar, P. et al. (2021). Severe Glutathione Deficiency, Oxidative Stress and Oxidant Damage in Adults Hospitalized with COVID-19: Implications for GlyNAC (Glycine and N-Acetylcysteine) Supplementation. *Antioxidants (Basel)*. **11**(1):50. doi: 10.3390/antiox11010050.



- De La Cruz Cortés, J.P. et al. (2021). Synergistic Effect of 3',4'-Dihidroxifenilglicol and Hydroxytyrosol on Oxidative and Nitrosative Stress and Some Cardiovascular Biomarkers in an Experimental Model of Type 1 Diabetes Mellitus. *Antioxidants*. **10**(12):1983. doi: 10.3390/antiox10121983.
- 14. Baxevanis, G.K. et al (2021). Tahini consumption improves metabolic and antioxidant status biomarkers in the postprandial state in healthy males. *Eur Food Res Technol*. doi: 10.1007/s00217-021-03828-5.
- Jacobson, M.H. et al. (2021). Organophosphate pesticides and progression of chronic kidney disease among children: A prospective cohort study. *Environ Int.* 155:106597. doi: 10.1016/j.envint.2021.106597.
- 16. Kumar, P. et al. (2021). Glycine and N-acetylcysteine (GlyNAC) supplementation in older adults improves glutathione deficiency, oxidative stress, mitochondrial dysfunction, inflammation, insulin resistance, endothelial dysfunction, genotoxicity, muscle strength, and cognition: Results of a pilot clinical trial. *Clin Transl Med.* **11**(3): e372. doi: 10.1002/ctm2.372.
- Szymańska, B. et al (2020). The Dependence between Urinary Levels of Angiogenesis Factors, 8-Iso-prostaglandin F2α, γ-Synuclein, and Interleukin-13 in Patients with Bladder Cancer: A Pilot Study. *J Oncol.* doi: 10.1155/2020/4848752.
- Headley, S.A. et al. (2020). The effects of 16-weeks of prebiotic supplementation and aerobic exercise training on inflammatory markers, oxidative stress, uremic toxins, and the microbiota in pre-dialysis kidney patients: a randomized controlled trial-protocol paper. *BMC Nephrol.* 21(1):517. doi: 10.1186/s12882-020-02177-x.
- 19. Jacobson, M.H. et al. (2020). Serially assessed bisphenol A and phthalate exposure and association with kidney function in children with chronic kidney disease in the US and Canada: A longitudinal cohort study. *PLoS Med.* **17**(10): e1003384. doi: 10.1371/journal.pmed.1003384.
- Agudelo, C.D. et al. (2020). Fermented Non-Digestible Fraction of Andean Berry (Vaccinium meridionale Swartz) Juice Induces Apoptosis in Colon Adenocarcinoma Cells. *Prev Nutr Food Sci.* 25(3):272-279. doi: 10.3746/pnf.2020.25.3.272.
- 21. Kumar, P. et al. (2020). Supplementing Glycine and N-acetylcysteine (GlyNAC) in Aging HIV Patients Improves Oxidative Stress, Mitochondrial Dysfunction, Inflammation, Endothelial Dysfunction, Insulin Resistance, Genotoxicity, Strength, and Cognition: Results of an Open-Label Clinical Trial. *Biomedicines*. 8(10): E390. doi: 10.3390/biomedicines8100390.
- 22. Gao, D. et al. (2020). In Vivo AAV Delivery of Glutathione Reductase Gene Attenuates Anti-aging Gene Klotho Deficiency-induced Kidney Damage. *Redox Biol.* doi: 10.1016/j.redox.2020.101692.
- 23. Marín-Echeverri, C. et al. (2020). Differential Effects of Agraz (Vaccinium meridionale Swartz) Consumption in Overweight and Obese Women with Metabolic Syndrome. *Journal of Food and Nutrition Research.* **8**(8):399-409. doi: 10.12691/jfnr-8-8-3.
- 24. Scarcello, E. et al. (2020). Amelioration of murine experimental colitis using biocompatible cyclosporine A lipid carriers. *Drug Deliv Transl Res.* doi: 10.1007/s13346-020-00835-z.
- 25. Wadsworth, D. et al. (2020). Randomised control study of oxidative stress in whole body vibration exercise. *JSES*. **4**(1):44-52. doi: 10.36905/jses.2020.01.07.
- Rawat, M. et al. (2020). Optimal Oxygen Targets in Term Lambs with Meconium Aspiration Syndrome and Pulmonary Hypertension. *Am J Respir Cell Mol Biol.* doi: 10.1165/rcmb.2019-0449OC.
- 27. Mistry, R.J. et al. (2020). Nicotinamide N-methyltransferase expression in SH-SY5Y human neuroblastoma cells decreases oxidative stress. *J Biochem Mol Toxicol*. doi: 10.1002/jbt.22439.



- 28. Rangarajan, S. et al. (2019). COX-2 derived prostaglandins as mediators of the deleterious effects of nicotine in chronic kidney disease. *Am J Physiol Renal Physiol*. doi: 10.1152/ajprenal.00407.2019.
- Ehnert-Russo, S.L. et al. (2019). Mercury Accumulation and Effects in the Brain of the Atlantic Sharpnose Shark (Rhizoprionodon terraenovae). *Arch Environ Contam Toxicol*. doi: 10.1007/s00244-019-00691-0.
- Sripetchwandee, J. et al. (2019). Deferiprone and efonidipine mitigated iron-overload induced neurotoxicity in wild-type and thalassemic mice. *Life Sci.* 239:116878. doi: 10.1016/j.lfs.2019.116878.

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: <u>tech@cellbiolabs.com</u> www.cellbiolabs.com

©2013-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.

