

FRAP[™] (Ferric Reducing Antioxidant Power) Detection Kit

Catalog No: K043-H1 (2 Plate)

Patent Protected

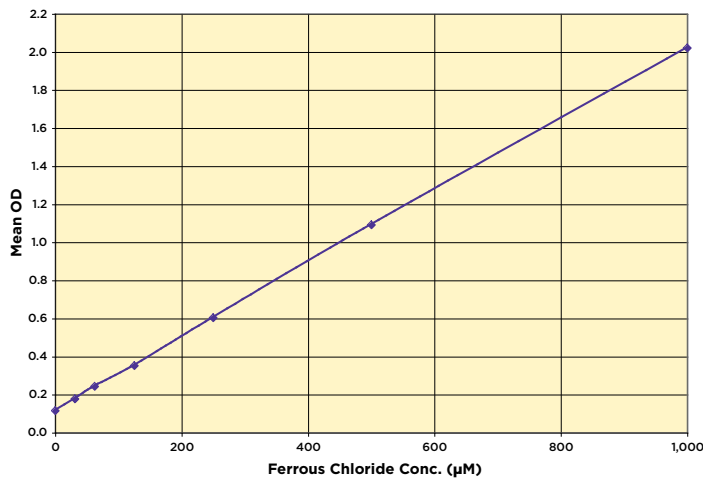
FEATURES

- ▶ Use Measure Ferric Reducing Anti-Oxidant Potential (FRAP)
- ▶ Sample Serum, Plasma, Urine, Food, Cosmetics, Additives
- ▶ Samples/Kit 89 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Time to Answer 30 Minutes
- ▶ Readout Colorimetric, 560 nm



SCIENTIFIC RELEVANCE

Potentially harmful reactive oxygen species (ROS) are produced as a consequence of normal aerobic metabolism. “Free Radicals” (FR) are usually removed or inactivated *in vivo* by a team of antioxidants. They are chemically stable atoms and molecules, which have one or more free electrons. Almost all biomolecules may be attacked by reactive free radicals. Free radicals are responsible for many pathological processes, or they can be generated as the result of the pathological stage and cause important secondary damage to biological systems and cells. Connections between free radicals and some serious diseases, including Parkinson’s and Alzheimer’s diseases, atherosclerosis, heart attacks, and chronic fatigue syndrome, have been demonstrated. However, short-term oxidative stress, the unbalance between the formation and scavenging of the reactive oxygen species, may be important in the prevention of aging due to triggering of the process known as mitohormesis. On average, 65 – 70% of the population is excessively impacted by oxidative stress caused by FRs.



Glutathione (GSH) Colorimetric Detection Kit

Catalog No: K006-H1 (4 Plate) K006-H1C-H/L (200 Cuvette)

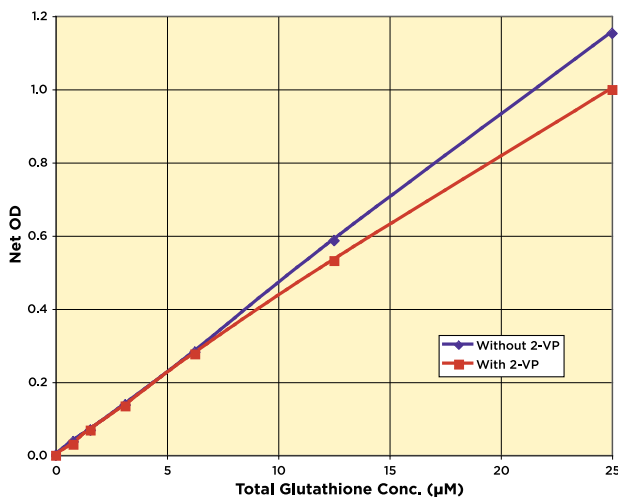
FEATURES

- ▶ Use Measure GSH/GSSG to Determine Oxidative Stress
- ▶ Sample Cells, RBCs, Serum, Plasma, Urine, and Tissue
- ▶ Sensitivity To 32 pmol/Sample
- ▶ Format 96-well or Cuvette
- ▶ Species Species Independent
- ▶ Samples/Kit 89 (Total and GSSG) in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 405 nm

**MULTI
SPECIES**

SCIENTIFIC RELEVANCE

Glutathione (L- γ -glutamyl-L-cysteinylglycine; GSH) is the highest concentration non-protein thiol in mammalian cells and is present in concentrations of 0.5 - 10 mM. GSH is a tripeptide that contains an unusual peptide linkage between the amine group of cysteine and the carboxyl group of the glutamate side-chain. It is an antioxidant, preventing damage to important cellular components caused by reactive oxygen species such as free radicals and peroxides. Glutathione reduces disulfide bonds formed within cytoplasmic proteins to cysteines by serving as an electron donor. In the process, glutathione is converted to its oxidized form, glutathione disulfide (GSSG). Glutathione is found mostly in its reduced form, since the enzyme that reverts it from its oxidized form, glutathione reductase, is constitutive and inducible upon oxidative stress. The ratio of reduced glutathione to oxidized glutathione within cells is often used as a measure of cellular toxicity.



Glutathione (GSH) Fluorescent Detection Kits

96 Well: **Catalog No: K006-F1 (1 Plate) K006-F5 (5 Plate)**
384 Well: **Catalog No: K006-F1D (2 Plate)**

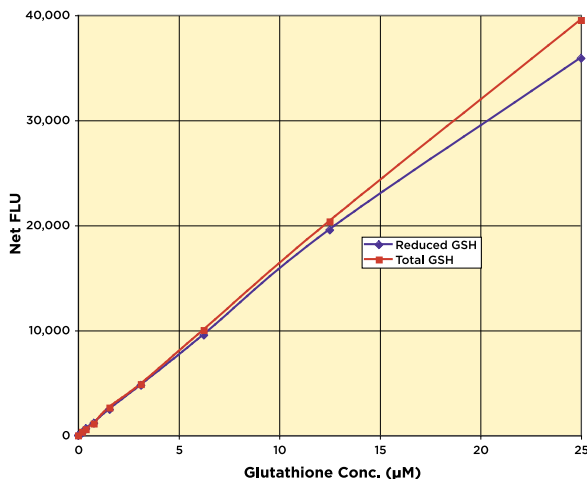
FEATURES

- ▶ **Use** Measure GSH/GSSG
- ▶ **Sample** Cells, RBC's, Serum, Plasma, Urine, and Tissues
- ▶ **Convenient** Measures Free and Total GSH in Same Sample Well
- ▶ **Species** Species Independent
- ▶ **Sensitivity** 45 nM Free GSH, 48 nM Total GSH
- ▶ **Samples/Kit** 96-well kits: 39 or 231 in Duplicate
 384-well kit: 183 in Duplicate
- ▶ **Stability** Liquid 4°C Stable Reagents
- ▶ **Readout** Fluorescent, 510 nm



SCIENTIFIC RELEVANCE

Glutathione (L-γ-glutamyl-L-cysteinylglycine; GSH) is the highest concentration non-protein thiol in mammalian cells and is present in concentrations of 0.5 - 10 mM. GSH is a tripeptide that contains an unusual peptide linkage between the amine group of cysteine and the carboxyl group of the glutamate side-chain. It is an antioxidant, preventing damage to important cellular components caused by reactive oxygen species such as free radicals and peroxides. Glutathione reduces disulfide bonds formed within cytoplasmic proteins to cysteines by serving as an electron donor. In the process, glutathione is converted to its oxidized form, glutathione disulfide (GSSG). Glutathione is found mostly in its reduced form, since the enzyme that reverts it from its oxidized form, glutathione reductase, is constitutive and inducible upon oxidative stress. The ratio of reduced glutathione to oxidized glutathione within cells is often used as a measure of cellular toxicity.



Glutathione Reductase Fluorescent Activity Kit

Catalog No: K009-F1 (1 Plate)

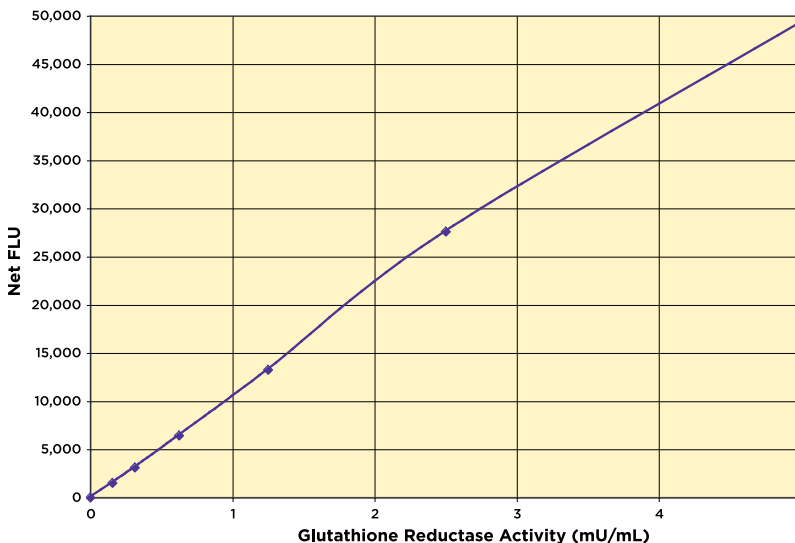
FEATURES

- ▶ Use Measure GR activity
- ▶ Sample RBCs, Serum, Plasma, and Cells
- ▶ Convenient 20 minute End Point or Kinetic Assay
- ▶ Sensitivity 9 µU/mL, World's Most Sensitive
- ▶ Species Species Independent
- ▶ Samples/Kit 41 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Fluorescent, 510 nm



SCIENTIFIC RELEVANCE

Glutathione reductase (GR) plays an indirect but essential role in the prevention of oxidative damage within the cell by helping to maintain appropriate levels of intracellular glutathione (GSH). GSH, in conjunction with the enzyme glutathione peroxidase (GP), is the acting reductant responsible for minimizing harmful hydrogen peroxide. The regeneration of GSH is catalyzed by GR. GR is a ubiquitous 100-120 kDa dimeric flavoprotein that catalyzes the reduction of oxidized glutathione (GSSG) to reduced glutathione, using β-nicotinamide dinucleotide phosphate (NADPH) as the hydrogen donor. NADPH has been suggested to also act as an indirectly operating antioxidant, given its role in the recycling of GSSG to GSH and thus maintaining the antioxidative power of glutathione.



Glutathione S-Transferase Fluorescent Activity Kit

Catalog No: K008-F1 (1 Plate)

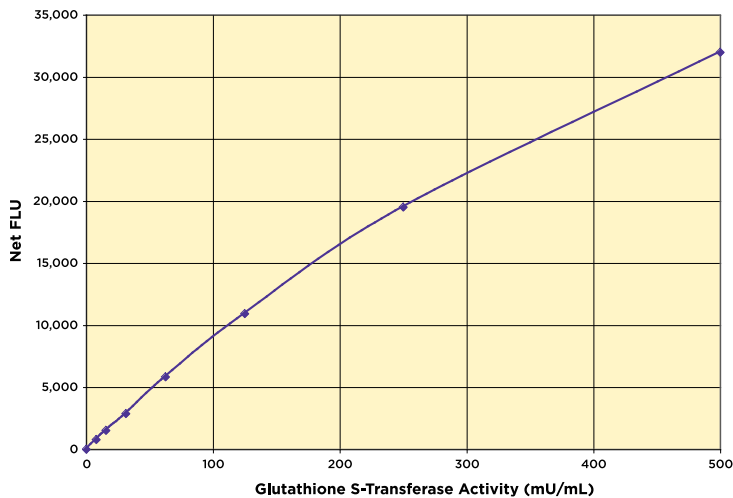
FEATURES

- ▶ Use Fluorescent Detection of GST Activity
- ▶ Sample Serum, Plasma, and Cell Lysates
- ▶ Samples/Kit 40 in Duplicate
- ▶ Convenient 30 Minute End Point or Kinetic Assay
- ▶ Sensitive < 100 μ U of GST Activity
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Fluorescent, 460 nm



SCIENTIFIC RELEVANCE

The Glutathione S-Transferase (GST) family of isozymes function to detoxify and neutralize a wide variety of electrophilic molecules by mediating their conjugation with reduced glutathione. Human GSTs are encoded by several gene families, and expressed in almost all tissues. Given its pivotal role in ameliorating oxidative stress/damage, GST activity has been repeatedly investigated as a biomarker for arthritis, asthma, COPD, and multiple forms of cancer, as well as an environmental marker. Examination of GST isoforms and activity in human cancers, tumors and tumor cell lines has revealed the predominance of the acidic pi class. Furthermore, this activity is thought to substantially contribute to the innate or acquired resistance of specific neoplasms to anticancer therapy.



Hydrogen Peroxide Colorimetric & Fluorescent Detection Kits

Colorimetric Catalog No: K034-H1 (2 Plate)

Fluorescent Catalog No: K034-F1 (2 Plate)

FEATURES

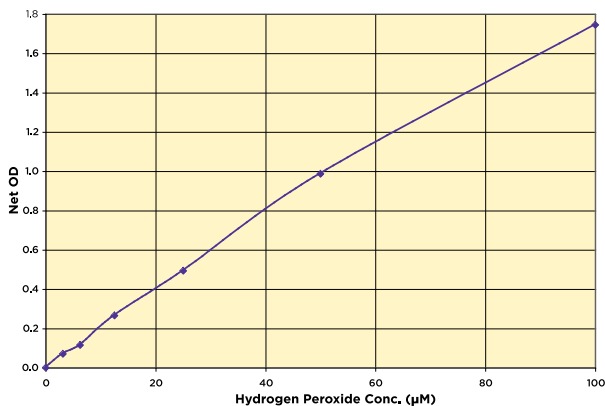
- ▶ Use Measure H_2O_2 in Any Sample
- ▶ Time to Answer 15 Minutes
- ▶ Sensitivity To 2 pmol (65 pg) H_2O_2
- ▶ Samples/Kit Colorimetric: 89 in Duplicate Fluorescent: 88 in Duplicate
- ▶ Readout Colorimetric: 560 nm Fluorescent: 585 nm



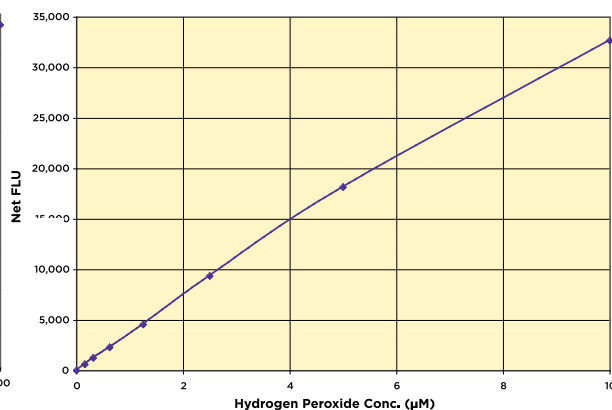
SCIENTIFIC RELEVANCE

In biological systems, incomplete reduction of O_2 during respiration produces superoxide anion ($O_2^{\cdot-}$), which is spontaneously or enzymatically dismutated by superoxide dismutase to H_2O_2 . Many cells produce low levels of $O_2^{\cdot-}$ and H_2O_2 in response to a variety of extracellular stimuli, such as cytokines (TGF- β 1, TNF- α , and various interleukins), peptide growth factors (PDGF; EGF, VEGF, bFGF, and insulin), the agonists of heterotrimeric G protein-coupled receptors (GPCR) such as angiotensin II, thrombin, lysophosphatidic acid, sphingosine 1-phosphate, histamine, and bradykinin, or by shear stress. The addition of exogenous H_2O_2 , or the intracellular production in response to receptor stimulation, affects the function of various proteins including protein kinases, protein phosphatases, transcription factors, phospholipases, ion channels, and G proteins. In 1894, Fenton described the oxidation of tartaric acid by Fe^{2+} and H_2O_2 , in which H_2O_2 and O_2 participate in the production of singlet oxygen and peroxyxynitrite. The generation of these species may be concurrent with reactions involving iron, and under some circumstances they might be important contributors to H_2O_2 toxicity.

Colorimetric Standard Curve



Fluorescent Standard Curve



Superoxide Dismutase (SOD) Colorimetric Activity Kit

Catalog No: K028-H1 (2 Plate)

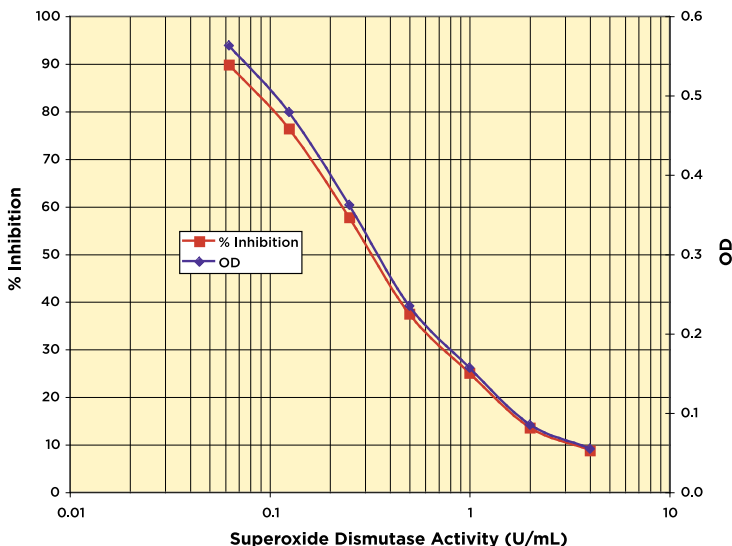
FEATURES

- ▶ Use Oxidative Stress Determination
- ▶ Sample Serum, Plasma, Urine, Cells and Tissue, and RBC
- ▶ Species Human and Other Mammalian Species
- ▶ Samples/Kit 89 in Duplicate
- ▶ Time to Answer 20 Minutes



SCIENTIFIC RELEVANCE

Short-lived and highly reactive oxygen species (ROS) such as $O_2^{\cdot -}$ (superoxide), $\cdot OH$ (hydroxyl radical), and H_2O_2 (hydrogen peroxide) are continuously generated *in vivo*. The cellular levels of ROS are controlled by antioxidant enzymes and small molecule antioxidants. The major antioxidant enzymes, superoxide dismutases (SODs), including copper-zinc superoxide dismutase (Cu/ZnSOD), manganese superoxide dismutase (MnSOD), and extracellular superoxide dismutase (EC-SOD). All play a critical roles in scavenging $O_2^{\cdot -}$. Decreased SOD activity results in elevated level of superoxide which in turn leads to decreased NO and increased peroxynitrite concentrations. The major intracellular SOD is a 32-kDa copper and zinc containing homodimer (Cu/Zn SOD). The mitochondrial SOD (MnSOD) is a manganese-containing 93-kDa homotetramer that is synthesized in the cytoplasm and translocated to the inner matrix of mitochondria. EC-SOD is the primary extracellular SOD enzyme and is highly expressed in many organs. Increased SOD activity levels are seen in Downs Syndrome, while decreased activity is seen in diabetes, Alzheimer’s disease, rheumatoid arthritis, Parkinson’s disease, uremic anemia, atherosclerosis, some cancers, and thyroid dysfunction.



Thiol Fluorescent Detection Kit

Catalog No: K005-F1 (1 Plate)

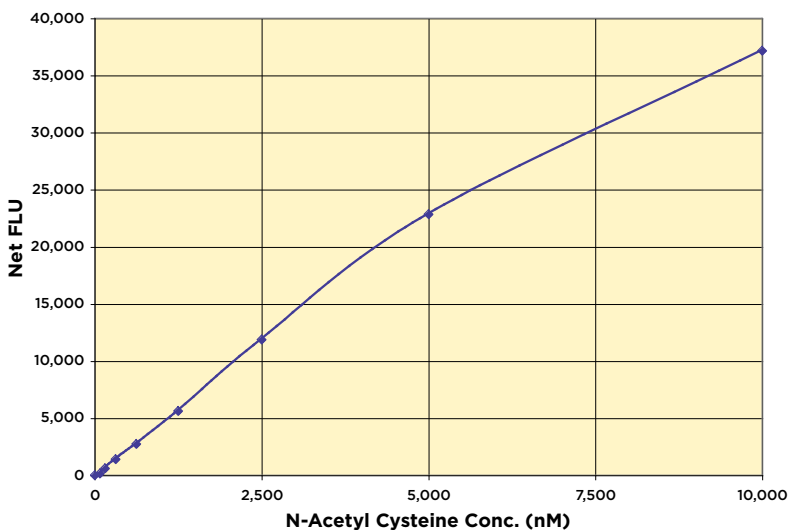
FEATURES

- ▶ Use Measure Thiol Content of Recombinant Proteins
- ▶ Adaptable Measure Protein Thiols (SH) in 6M GuHCl Buffers
- ▶ Sensitivity 4.62 nM
- ▶ Time to Answer 30 Minutes
- ▶ Samples/Kit 39 in Duplicate
- ▶ Stability Non-Toxic, Liquid 4°C Stable Reagents
- ▶ Readout Fluorescent, 510 nm



SCIENTIFIC RELEVANCE

Free thiols in biological systems have important roles. Oxidatively modified thiol groups of cysteine residues are known to modulate the activity of a growing number of proteins. One of the most pressing problems is to accurately determine the extent of modification of specific amino acids, such as cysteine residues, in a complex protein sample, especially in the presence of chaotropic agents such as guanidine hydrochloride. Typical methods such as using Ellman's reagent have limited sensitivity requiring large quantities of purified recombinant or native protein.





www.ArborAssays.com
(734) 677-1774
info@ArborAssays.com
©2014

180403

WEB 180403