



ARBOR ASSAYS
Interactive Assay Solutions™



Table of Contents

DetectX® Assay Kits	Page
Adrenocorticotrophic Hormone (ACTH) ELISA Kits	4
Catalase Colorimetric and Fluorescent Activity Kits	5
Ceruloplasmin Colorimetric Activity Kit	6
Corticosterone Chemiluminescent ELISA Kits	7
Corticosterone Multi-Format ELISA Kits	8
Cortisol ELISA Kits	9
Cortisone Chemiluminescent ELISA Kits	10
Cortisone ELISA Kits	11
FRAP™ (Ferric Reducing Antioxidant Power) Detection Kit	12
Glutathione Colorimetric Detection Kits	13
Glutathione Fluorescent Detection Kits	14
Glutathione Reductase Fluorescent Activity Kit	15
Glutathione S-Transferase (GST) Fluorescent Activity Kit	16
Hydrogen Peroxide (H ₂ O ₂) Colorimetric and Fluorescent Detection Kits	17
Superoxide Dismutase (SOD) Colorimetric Activity Kit	18
Thiol Fluorescent Detection Kit	19



ONLINE:

www.ArborAssays.com/order-form

PHONE:

Call 734-677-1774 or Toll Free: 855-677-1774. Monday-Friday 8:30 am to 5:30 pm, EST.

FAX:

Send faxes to 734-677-6860.

E-MAIL:

Send E-mail orders to Orders@ArborAssays.com

DISTRIBUTORS:

Check our website at www.ArborAssays.com/distributors for a list of distributors.

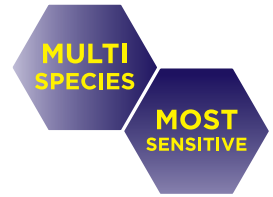


Adrenocorticotrophic Hormone (ACTH) ELISA Kits

K072-H1 (1 Plate) | K072-H5 (5 Plate)

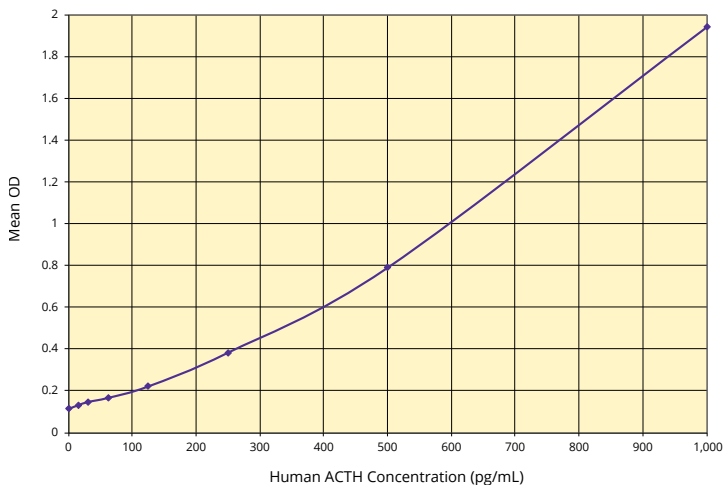
FEATURES

- ▶ Use Measure ACTH in Plasma Samples
- ▶ Sample Plasma
- ▶ Sensitivity 12.71 pg/mL
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Adrenocorticotrophic hormone (ACTH, also adrenocorticotropin, corticotropin) is an important component of the hypothalamic-pituitary-adrenal axis (HPA) and is produced in response to biological stress. Its principal effects are increased production and release of glucocorticoids (GCs). Stress-induced secretion of the peptide hormone Corticotropin Releasing Hormone (CRH) stimulates pituitary ACTH secretion. Circulating ACTH binds to melanocortin receptors on the surface of adrenal zona cells, inducing the synthesis and release of all adrenal steroids, aldosterone, GCs and adrenal androgens. ACTH is also the principal modulator of cortisol and corticosterone. In addition to the stress response, ACTH synthesis is related to the circadian rhythm in many organisms. Measurement of plasma ACTH is helpful in the differential diagnosis of pituitary Cushing's disease, Addison's disease, adrenal tumors, adrenal hyperplasia, and ectopic ACTH syndrome.



Catalase Colorimetric Activity Kits

Colorimetric: K033-H1 (2 Plate)
 Fluorescent: K033-F1 (2 Plate)

FEATURES

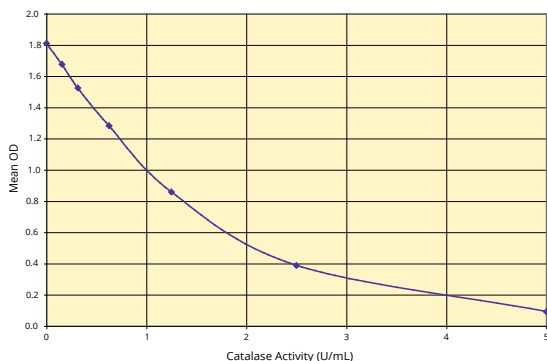
- ▶ Use Measure Catalase Activity in Any Sample
- ▶ Time to Answer 45 Minutes
- ▶ Sensitivity Measure as Little as 0.052 U/mL
- ▶ Samples/Kit 89 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Species Species Independent
- ▶ Readout Colorimetric, 560 nm
 Fluorescent, 590 nm em/520 nm ex



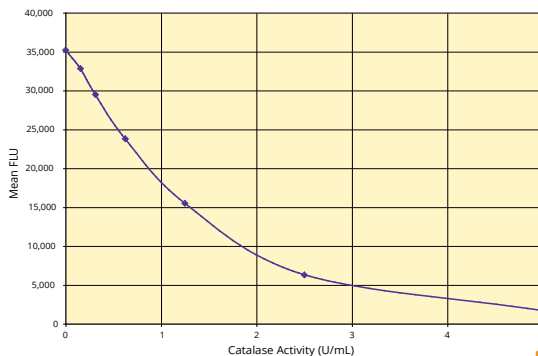
SCIENTIFIC RELEVANCE

Hydrogen peroxide, (H₂O₂) is one of the most frequently occurring reactive oxygen species. It is formed either in the environment, as a by-product of aerobic metabolism, superoxide formation and dismutation, or as a product of oxidase activity. Both excessive hydrogen peroxide and its decomposition product hydroxyl radical, are harmful for most cell components. Its rapid removal is essential for all aerobically living prokaryotic and eukaryotic cells. One of the most efficient ways of removing peroxide is through the enzyme catalase, which is encoded by a single gene and is highly conserved among species. Mammals, including humans and mice, express catalase in all tissues. A high concentration of catalase can be found in the liver, kidneys and erythrocytes. The expression is regulated at transcription, post-transcription and post-translation levels. High catalase activity is detected in peroxisomes.

Colorimetric



Fluorescent



Ceruloplasmin (Cp) Colorimetric Activity Kit

K035-H1 (2 Plate)

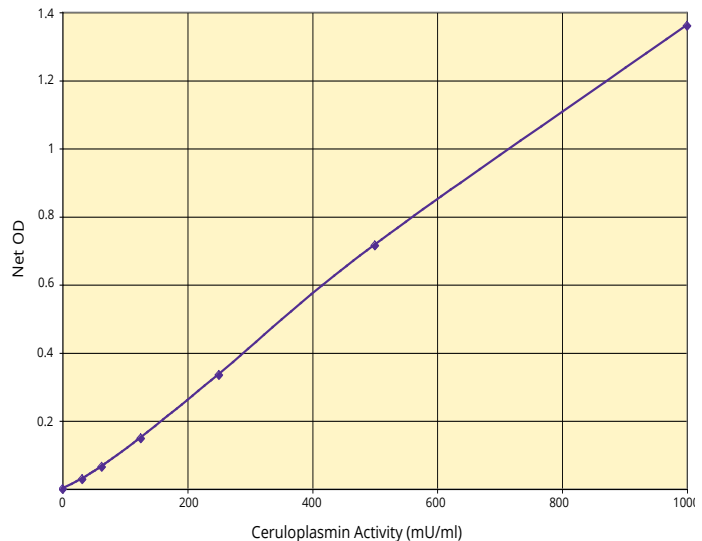
FEATURES

- ▶ Use Non-Invasive Pregnancy Marker
- ▶ Sample Urine, Serum
- ▶ Validation Humans, Felids, Polar Bear, Panda
- ▶ Species Multiple Species
- ▶ Time to Answer 60 Minutes
- ▶ Format 96-Well
- ▶ Samples/Kit 89 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 560 nm



SCIENTIFIC RELEVANCE

Ceruloplasmin (Cp) is an acute phase multicopper oxidase enzyme that normally plays a protective role in responses to immune-provoking stimuli and is also associated with reproduction. Estrogens alter the subcellular distribution of copper in the liver, leading to an increase in plasma copper levels and subsequent ceruloplasmin synthesis. Serum levels of Cp have been shown to increase during normal pregnancy in some species possibly as a protection against the oxidative costs of reproduction. In giant pandas and some felids, urinary Cp activity has been shown to be elevated in pregnant vs. pseudopregnant animals beginning in the first week of gestation and continuing throughout the luteal phase.



Corticosterone Chemiluminescent ELISA Kits

K014-C1 (1 Plate) | K014-C5 (5 Plate)

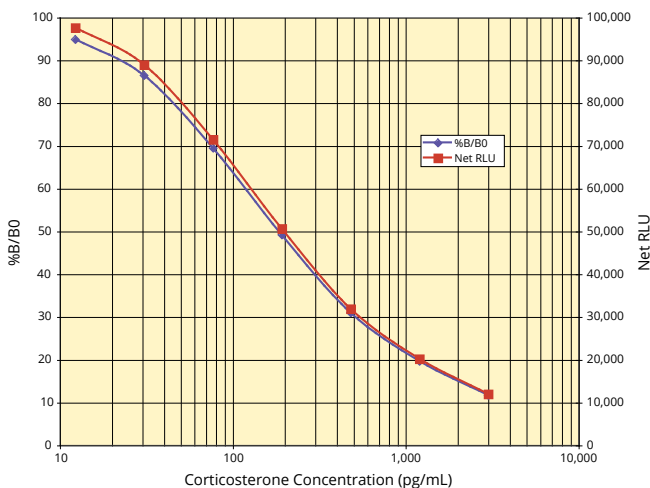
FEATURES

- ▶ Use Stress Marker in < 2µL Serum or Plasma
- ▶ Sample Serum, Plasma, Urine, Respiratory Vapor, TCM. and extracts of Feces, Feathers and Hair
- ▶ Validation Mice, Rats, Humans, Primates, Birds, Cats, Ungulates, Whale, Lizard
- ▶ Sensitivity 6.7 pg/mL
- ▶ Time to Answer 2 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 39 or 231 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Glow Luminescent



SCIENTIFIC RELEVANCE

Corticosterone (Kendall's Compound 'B') is a glucocorticoid secreted by the cortex of the adrenal gland. It is produced in response to stimulation of the adrenal cortex by ACTH and is the precursor of aldosterone. Corticosterone is a major indicator of stress and is the major stress steroid produced in non-human mammals. Studies involving corticosterone and levels of stress have found evidence of impairment of long term memory retrieval, chronic corticosterone elevation due to dietary restrictions, and corticosterone elevation in response to burn injuries. In addition to stress levels, corticosterone is believed to play a decisive role in sleep-wake patterns.



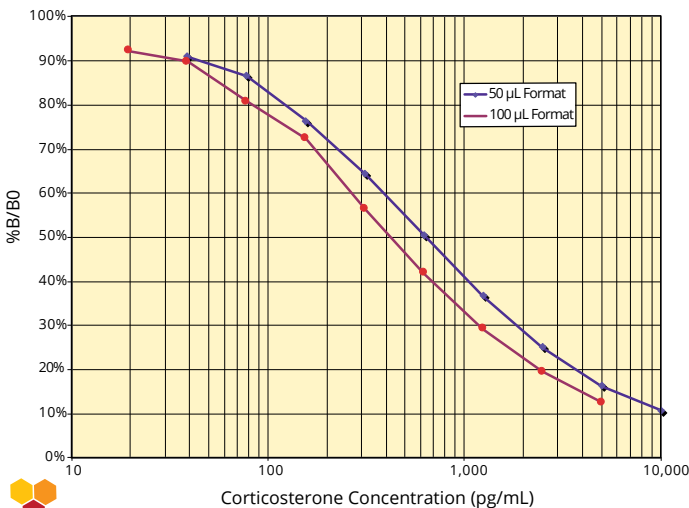
Corticosterone Multi-Format ELISA Kits

K014-H1 (1 Plate) | K014-H5 (5 Plate)

K014-H1W (1 Whole Plate) | K014-H5W (5 Whole Plate)

FEATURES

- ▶ Use Stress Marker in < 2 μ L Serum or Plasma
- ▶ Sample Serum, Plasma, Urine, Respiratory Vapor, TCM and Feather, Hair and Fecal Extracts
- ▶ Multi-Format Standard range of 10,000–39.06 pg/mL or 5,000–19.53 pg/mL
- ▶ Sensitivity 50 μ L Format: 20.9 pg/mL
100 μ L Format: 17.5 pg/mL
- ▶ Validation Mice, Rats, Humans, Primates, Birds, Cats, Ungulates, Whale, Lizard
- ▶ Time to Answer 1.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 37 or 229 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Corticosterone (Kendall's Compound 'B') is a glucocorticoid secreted by the cortex of the adrenal gland. It is produced in response to stimulation of the adrenal cortex by ACTH and is the precursor of aldosterone. Corticosterone is a major indicator of stress and is the major stress steroid produced in non-human mammals. Studies involving corticosterone and levels of stress have found evidence of impairment of long term memory retrieval, chronic corticosterone elevation due to dietary restrictions, and corticosterone elevation in response to burn injuries. In addition to stress levels, corticosterone is believed to play a decisive role in sleep-wake patterns.



Cortisol ELISA Kits

K003-H1 (1 Strip Plate) | K003-H5 (5 Strip Plate)
 K003-H1W (1 Whole Plate) | K003-H5W (5 Whole Plate)

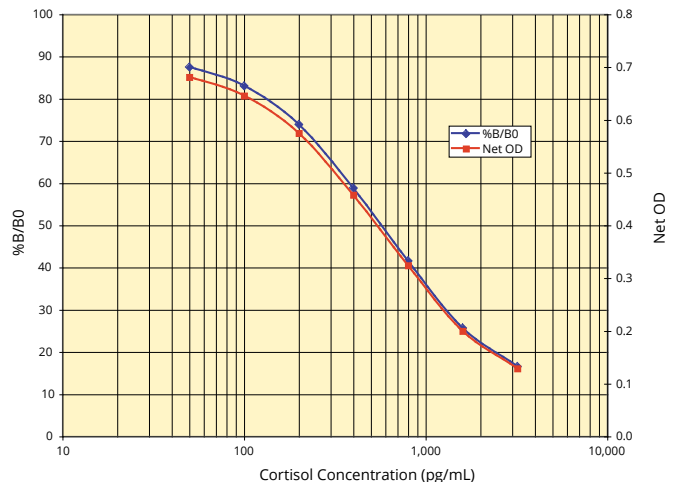
FEATURES

- ▶ Use Stress Marker in as Little as 1 μ L Serum or Plasma
- ▶ Sample Serum, Plasma, Saliva, Urine, Respiratory Vapor, and Hair and Fecal Extracts
- ▶ Validation Rodent, Primates, Ungulates, Fish, Whale, Canine
- ▶ Time to Answer 1.5 Hours
- ▶ Format 96-Well, Break-Apart Strip or Whole Plates
- ▶ Species Species Independent
- ▶ Samples/Kit 39 or 231 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Cortisol (hydrocortisone, Kendall’s Compound ‘F’) is the primary glucocorticoid produced and secreted by the adrenal cortex. It is often referred to as the “stress hormone” as it affects blood pressure, blood sugar levels, and other actions of stress adaptation. Immunologically, cortisol functions as an important anti-inflammatory and plays a role in hypersensitivity, immunosuppression, and disease resistance. In the metabolic aspect, cortisol promotes gluconeogenesis, liver glycogen deposition, and the reduction of glucose utilization. Production of cortisol follows an ACTH-dependent circadian rhythm, with a peak level in the morning and decreasing levels throughout the day. All but 4% of serum cortisol is bound to proteins including corticosteroid binding globulin and serum albumin. Abnormal cortisol levels are being evaluated for correlation with a variety of different conditions, such as prostate cancer, depression, schizophrenia, Cushing’s Syndrome, and Addison’s Disease.

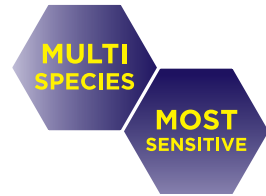


Cortisone Chemiluminescent ELISA Kits

K017-C1 (1 Plate) | K017-C5 (5 Plate)

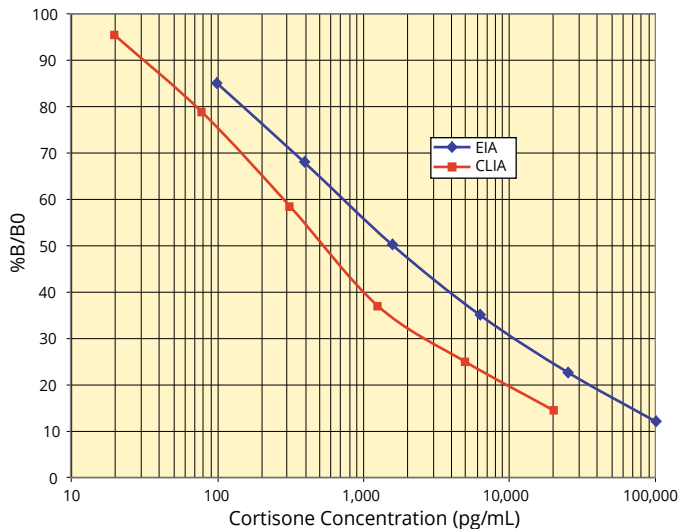
FEATURES

- ▶ Use Stress Marker in as Little as 1 μ L Serum or Plasma
- ▶ Sample Serum, Plasma, Urine, Saliva, Fecal Extracts
- ▶ Validation Mice, Rats, Humans, Monkeys, Felids, Ungulates, Canine
- ▶ Time to Answer 2 Hours
- ▶ Species Species Independent
- ▶ Samples/Kit 37 or 229 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Glow Luminescent



SCIENTIFIC RELEVANCE

Cortisone (C₂₁H₂₈O₅, Kendall's Compound 'E') was identified by extraction from bovine suprarenal gland tissue. Cortisol and cortisone concentrations vary due to the activity of two 11 β -hydroxysteroid dehydrogenases (11 β -HSD). 11 β -HSD1 is found primarily in the liver where it converts cortisone to cortisol while 11 β -HSD2 is found in tissues such as the kidney where cortisol receptor binding is required. This glucocorticoid "shuttle" helps to initiate and regulate the anti-inflammatory response.



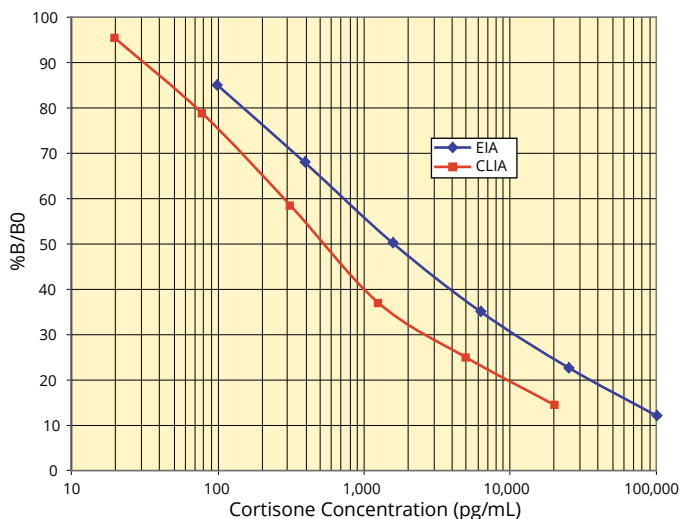
FEATURES

- ▶ Use Stress Marker in as Little as 1 µL Serum or Plasma
- ▶ Sample Serum, Plasma, Urine, Saliva, Fecal Extracts
- ▶ Validation Mice, Rats, Humans, Monkeys, Felids, Ungulates, Canine
- ▶ Time to Answer 2.5 Hours
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Cortisone (C₂₁H₂₈O₅, Kendall’s Compound ‘E’) was identified by extraction from bovine suprarenal gland tissue. Cortisol and cortisone concentrations vary due to the activity of two 11β-hydroxysteroid dehydrogenases (11β-HSD). 11β-HSD1 is found primarily in the liver where it converts cortisone to cortisol while 11β-HSD2 is found in tissues such as the kidney where cortisol receptor binding is required. This glucocorticoid “shuttle” helps to initiate and regulate the anti-inflammatory response.



FRAP™ (Ferric Reducing Antioxidant Power) Detection Kit

K043-H1 (2 Plate)

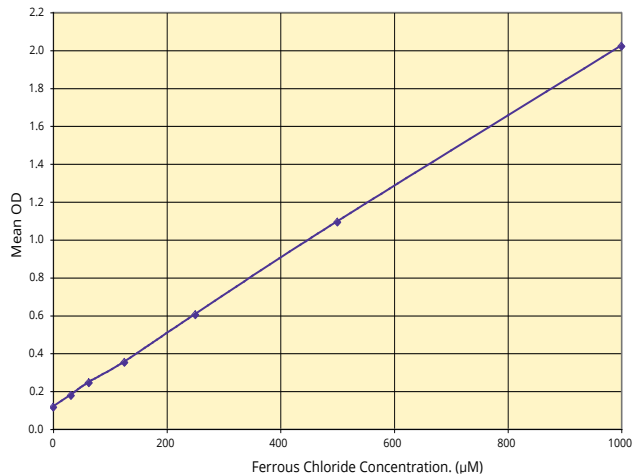
FEATURES

- ▶ Use Measure Ferric Reducing Anti-Oxidant Potential (FRAP) of Samples
- ▶ Samples Serum, Plasma, Tissue, Saliva, Cell Lysates, 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 preventing 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 Kits

K006-H1 (4 Plate)

K006-H1C-H/L (200 Cuvette)

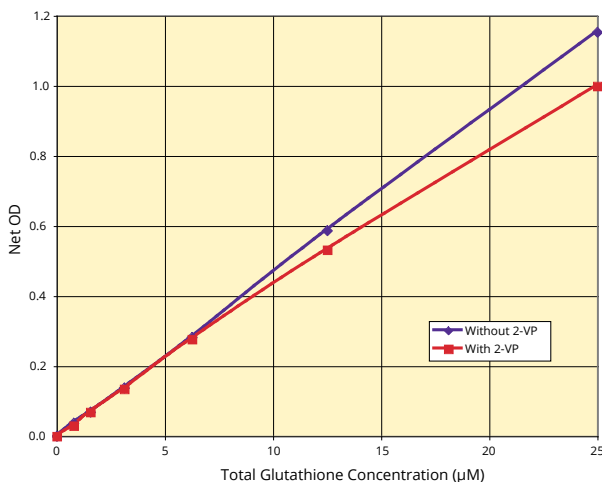
FEATURES

- ▶ Use Measure Total GSH and GSSG to Determine Oxidative Stress
- ▶ Samples Whole Blood, Serum, Plasma, Erythrocytes, Urine, Lysates, TCM
- ▶ Sensitivity 0.634 μ M (Plate-based Format)
- ▶ Format 96-Well or Cuvette
- ▶ Species Species Independent
- ▶ Samples/Kit K006-H1: 89 (Total and GSSG) in Duplicate
K006-H1C-H/L: 43 (Total and GSSG) in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 405 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 (GSH) Fluorescent Detection Kits

K006-F1 (1 Plate) | K006-F5 (5 Plate)
 K006-F1D (384-Well Plate)

FEATURES

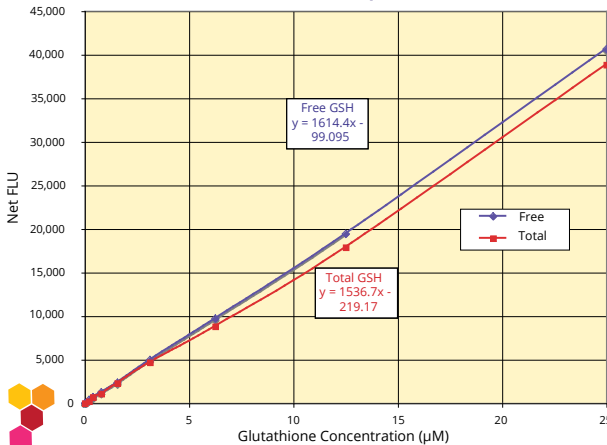
- ▶ Use: Measure GSH/GSSG to Determine Oxidative Stress
- ▶ Sample: Whole Blood, Serum, Plasma, Erythrocytes, Urine, Lysates, TCM
- ▶ Species: Species Independent
- ▶ Sensitivity: 45 nM Free GSH, 48 nM Total GSH
- ▶ Samples/Kit: 39 or 231 in Duplicate (K006-F1/F5)
183 in Duplicate (K006-F1D)
- ▶ Stability: Liquid 4°C Stable Reagents
- ▶ Readout: Fluorescent, 510 nm em/370-410 nm ex



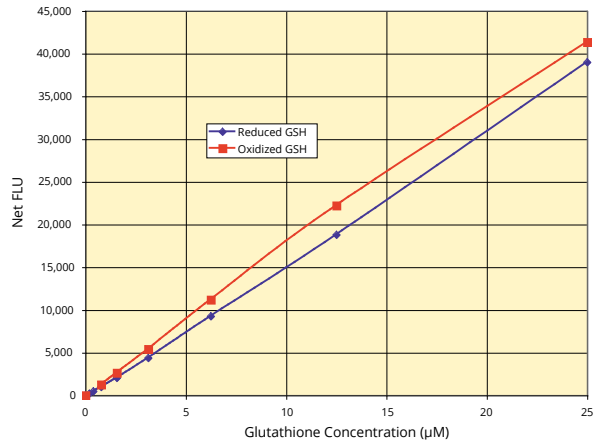
SCIENTIFIC RELEVANCE

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.

K006-F1/F5



K006-F1D



Glutathione Reductase Fluorescent Activity Kit

K009-F1 (1 Plate)

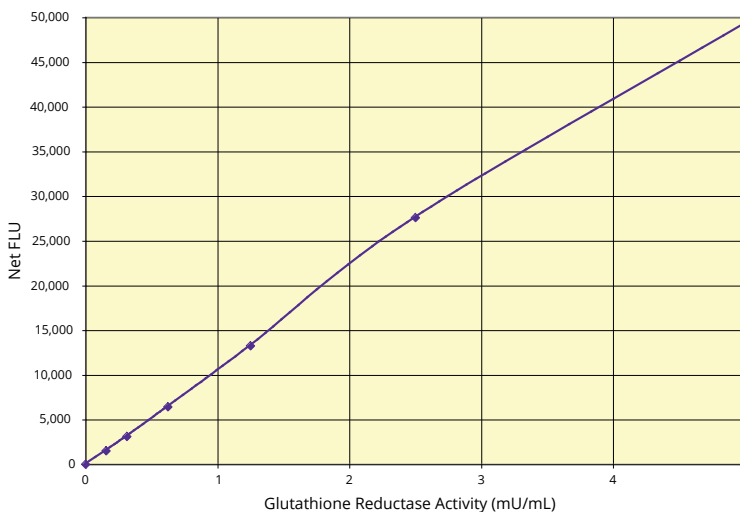
FEATURES

- ▶ Use Measure GR activity
- ▶ Sample RBCs, Serum, Plasma, Cells, Tissues
- ▶ Convenient 20 minute End Point or Kinetic Assay
- ▶ Sensitivity 9 μ U/mL, World's Most Sensitive
- ▶ Species Species Independent
- ▶ Samples/Kit 41 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Fluorescent, 510 nm em/370-410 nm ex



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

K008-F1 (1 Plate)

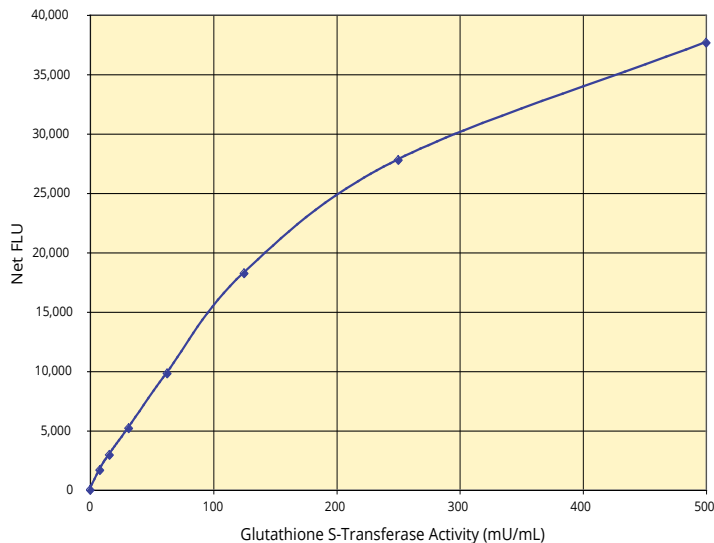
FEATURES

- ▶ Use Measure GST Activity
- ▶ Sample Serum, Plasma, Urine, Cell Lysates
- ▶ Samples/Kit 40 in Duplicate
- ▶ Convenient 30 Minute End Point or Kinetic Assay
- ▶ Sensitivity < 100 μ U of GST Activity
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Fluorescent, 460 nm em/370-410 nm ex



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 five gene families, expressing in almost all tissues as four cytosolic and one microsomal forms. 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: K034-H1 (2 Plate)

Fluorescent: K034-F1 (2 Plate)

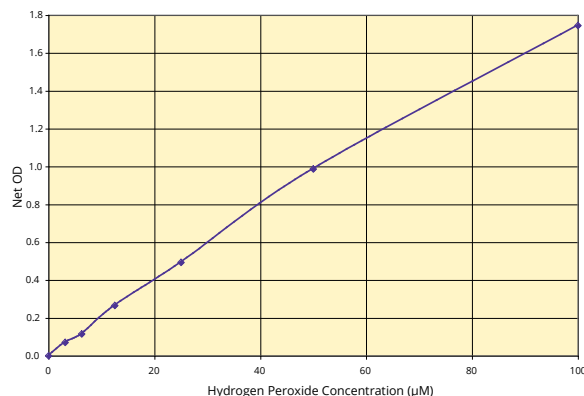
FEATURES

- ▶ Use Measure H_2O_2 in Any Sample
- ▶ Sample Urine, Buffer, TCM
- ▶ Rapid 15 Minutes
- ▶ Sensitive Colorimetric: 91.3 pmol (310 ng)
Fluorescent: < 2 pmole (65 pg)
- ▶ Samples/Kit Colorimetric: 89 in Duplicate
Fluorescent: 88 in Duplicate
- ▶ Readout Colorimetric: 560 nm
Fluorescent: 590 nm em/520 nm ex

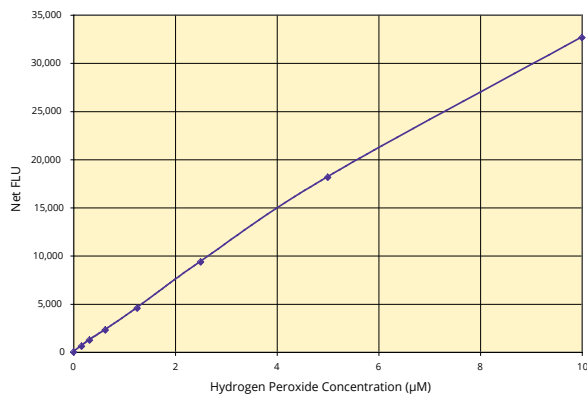
SCIENTIFIC RELEVANCE

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, and 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 . H_2O_2 and O_2 may participate in the production of singlet oxygen and peroxyxynitrite and the generation of these species may be concurrent with reactions involving iron, which under some circumstances might be important contributors to H_2O_2 toxicity.

Colorimetric



Fluorescent



Superoxide Dismutase (SOD) Activity Kit

K028-H1 (2 Plate)

FEATURES

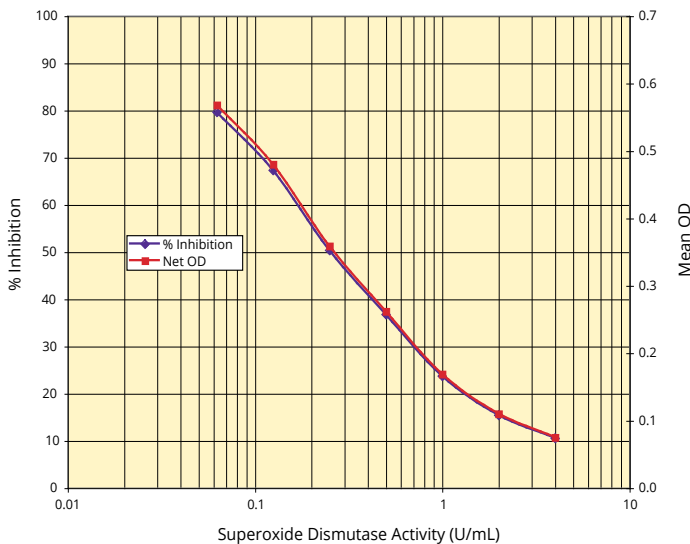
- ▶ Use Oxidative Stress Determination
- ▶ Sample Serum, Plasma, Cells, Tissue Buffers, Erythrocytes
- ▶ Species Human and Other Mammalian Species
- ▶ Samples/Kit 88 in Duplicate
- ▶ Time to Answer 20 Minutes
- ▶ Readout Colorimetric, 450 nm



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

K005-F1 (1 Plate)

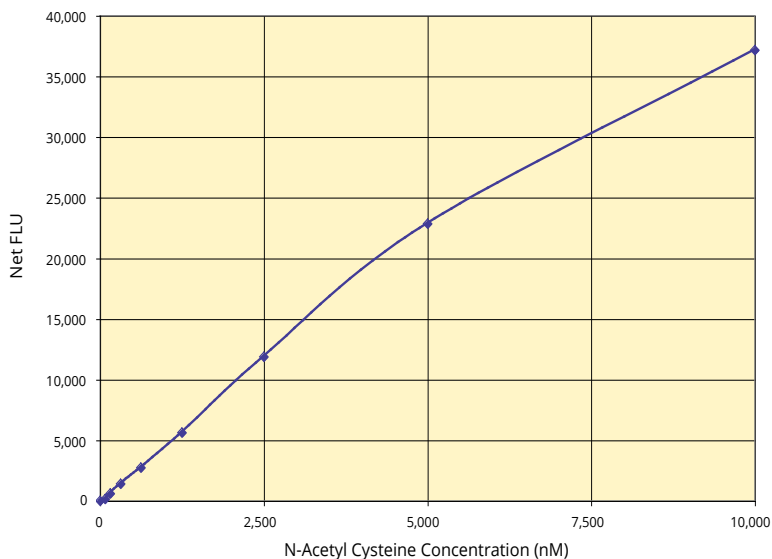
FEATURES

- ▶ Use Measure Thiol Content of Proteins and Peptides
- ▶ Adaptable Measure SH easily in 8M GuHCl Buffers
- ▶ Sensitivity 4.62 nM
- ▶ Time to Answer 30 Minutes
- ▶ Species Species Independent
- ▶ Samples/Kit 39 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Fluorescent, 510 nm em/370-410 nm ex



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. As such, it is important to be able to accurately determine the extent of modification of specific amino acids, such as cysteine residues. This is especially difficult in a complex protein sample, especially in the presence of chaotropic agents such as guanidine hydrochloride. Typical methods using Ellman's reagent do not have sufficient sensitivity to allow economical detection of free SH groups.





STRESS

REPRODUCTIVE

OXIDATIVE
STRESS



METABOLISM

VOLUME &
NORMALIZATION



KIDNEY
INJURY



INFLAMMATION



CELL
SIGNALING



ARBOR
ASSAYS™