



ARBOR ASSAYS
Interactive Assay Solutions™

REPRODUCTIVE

Table of Contents

DetectX® Assay Kits	Page
17-Hydroxyprogesterone ELISA Kits	4
Aldosterone ELISA & Chemiluminescent ELISA Kits	5
Allopregnanolone ELISA Kits	6
Androstenedione ELISA Kits	7
Arg8-Vasopressin (AVP) ELISA & Chemiluminescent ELISA Kits	8
Ceruloplasmin (Cp) Colorimetric Activity Kit	9
Creatinine Urinary Detection Kits/Solution	10
Dehydroepiandrosterone Sulfate (DHEA-S) ELISA Kits	11
Epiandrosterone ELISA Kits	12
Estradiol ELISA Kits	13
Serum Estradiol ELISA Kits	14
Estriol ELISA Kits	15
Estrone ELISA Kits	16
Estrone-3-Glucuronide (E1G) ELISA Kits	17
Estrone-3-Sulfate (E1S) ELISA Kits	18
Levonorgestrel (LNG) ELISA Kits	19
Oxytocin ELISA & Chemiluminescent ELISA Kits/Solutions	20
PGFM (13,14-dihydro-15-keto-Prostaglandin F2a) ELISA Kits	21
Pregnanediol-3-Glucuronide (PDG) ELISA Kits	22
Progesterone ELISA Kits	23
Progesterone Metabolites ELISA Kits	24
Prolactin (PRL) ELISA Kit	25
Testosterone ELISA Kits	26
Testosterone ELISA Kits - Improved Sensitivity	27



ONLINE:

www.ArborAssays.com/order-form

PHONE:

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

FAX:

Send faxes to [734-677-6860](tel: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.



17-Hydroxyprogesterone ELISA Kits

K053-H1 (1 Plate) | K053-H5 (5 Plate)

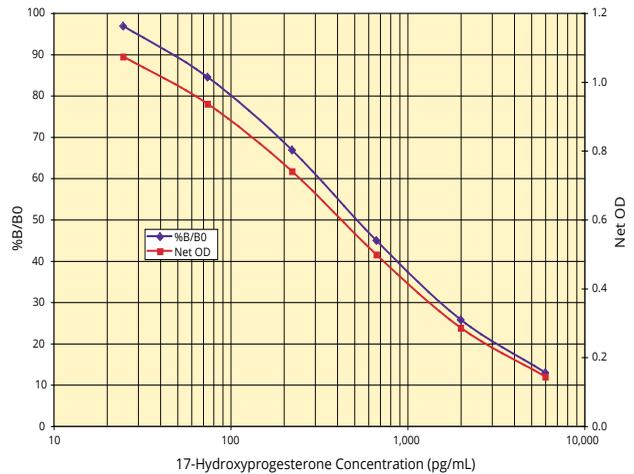
FEATURES

- ▶ Use Congenital Adrenal Hyperplasia Marker
- ▶ Sample Urine and Extracted Serum, Plasma, and Fecal
- ▶ Time to Answer 1.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

17-Hydroxyprogesterone is a steroid hormone belonging to the androgen group and is found and is in mammals, reptiles, birds, and other vertebrates. It was first isolated by Pfiffner and North in 1940. It is primarily produced in the adrenal glands but is also produced in the corpus luteum of the ovary. It is hydroxylated at the 11 and 21 positions to produce cortisol. A deficiency of either 11- or 21-hydroxylase results in decreased cortisol synthesis, and feedback inhibition of adrenocorticotrophic hormone (ACTH) secretion is lost. Consequently increased pituitary release of ACTH will increase production of 17HO-P. However, if 17-alpha-hydroxylase or 3β-hydroxysteroid dehydrogenase type 2 is deficient, then 17HO-P levels will be low with either a possible increase in progesterone or pregnenolone respectively. Normal levels are 3-90 ng/dL in children. In women, normal levels are 20-100 ng/dL prior to ovulation, and 100-500 ng/dL during the luteal phase.



Aldosterone ELISA & Chemiluminescent ELISA Kits

ELISA: K052-H1 (1 Plate) | K052-H5 (5 Plate)

Chemiluminescent ELISA: K052-C1 (1 Plate) | K052-C5 (5 Plate)

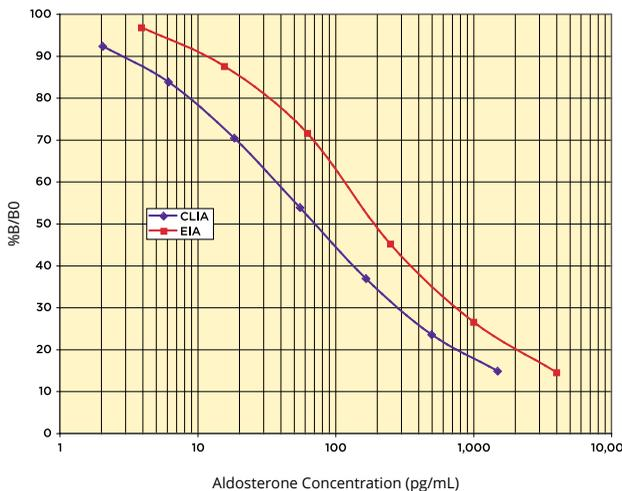
FEATURES

- ▶ Use Reproductive Assessment
- ▶ Sample Urine and Extracted Serum, Plasma, and Dried Feces
- ▶ Time to Answer ELISA: 2.5 Hours
Chemiluminescent: Overnight
- ▶ Sensitivity ELISA: 4.97 pg/mL
Chemiluminescent: 1.84 pg/mL
- ▶ Samples/Kit ELISA: 40 or 232 in Duplicate
Chemiluminescent: 39 or 231 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout ELISA: Colorimetric, 450 nm
Chemiluminescent: Glow Luminescent



SCIENTIFIC RELEVANCE

Aldosterone is a mineralocorticoid first isolated by the husband and wife team of Simpson and Tait at University College, London. Aldosterone controls the sodium-potassium balance through unidirectional salt reabsorption in a variety of tissues and glands. Synthesized from cholesterol in the zona glomerulosa of the adrenal cortex, secretion is regulated through the renin-angiotensin system. Angiotensin II and potassium stimulate primary secretion by increasing the rate of production of the steroid. Peripheral aldosterone levels are dependent on age and body position. Aldosterone measurement is useful in the investigation of primary and secondary aldosteronism including vascular disease, salt depletion, potassium loading, cardiac failure with ascites, and pregnancy.



Allopregnanolone ELISA Kits

K061-H1 (1 Plate) | K061-H5 (5 Plate)

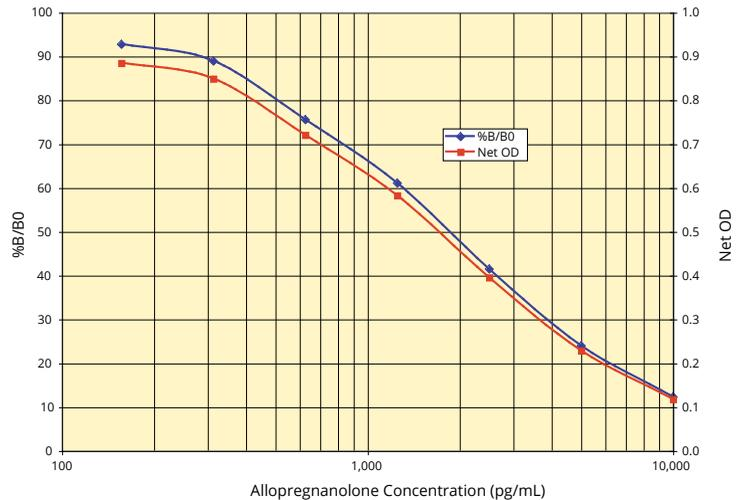
FEATURES

- ▶ Use Measure allopregnanolone in a variety of samples
- ▶ Sample Urine, TCM or Extracted Serum, Plasma, Dried Feces
- ▶ Time to Answer 2.5 Hours or Overnight
- ▶ Range 10,000 – 156.3 pg/mL
- ▶ Species Species Independent
- ▶ Samples/Kit 39 or 231 in Duplicate
- ▶ Cross Reactivity Low Cross Reactivity to Progesterone and Metabolites
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Allopregnanolone (3 α -hydroxy-5 α -pregnan-20-one, THP, THPROG) is a prototypic neurosteroid present in the blood and the brain. It is a metabolite of progesterone and potent modulator of GABA_A receptors. Allopregnanolone has pharmacological properties including anxiolytic and anticonvulsant activity. The biosynthesis of allopregnanolone involves the conversion of progesterone into 5 α -dihydroprogesterone by the enzyme 5 α -reductase type I. Subsequently, 3 α -hydroxysteroid oxidoreductase isoenzymes convert this intermediate into allopregnanolone. Anxiety and depression are common side effects of 5 α -reductase inhibitors such as finasteride and dutasteride, and they are believed to be caused, in part, by the prevention of the endogenous production of allopregnanolone. Allopregnanolone aids neurogenesis and has been found to reverse neuron proliferative deficit and cognitive deficits in mouse models of Alzheimer's disease.



FEATURES

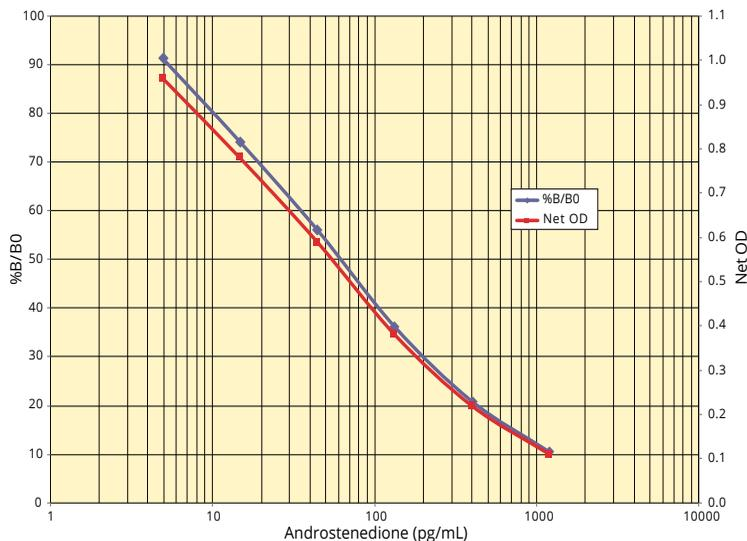
- ▶ Use Assessment of sexual development disruption; marker of androgen biosynthesis
- ▶ Sample Saliva, Urine and Extracted Serum or Plasma
- ▶ Time to Answer 2.5 Hours
- ▶ Sensitivity 2.3 pg/mL
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



MULTI
SPECIES

SCIENTIFIC RELEVANCE

Androstenedione is a steroid hormone also known as androst-4-ene-3,17-dione. It is synthesized from dehydroepiandrosterone (DHEA) or 17-hydroxyprogesterone. By itself, androstenedione is a weak androgen though it is a precursor for biosynthesis of stronger androgens, such as testosterone and estrogens. Therefore, androstenedione is classified as a controlled drug and a supplement to enhance athletic performance and body energy.



Arg⁸-Vasopressin (AVP) ELISA & Chemiluminescent ELISA Kits

ELISA: K049-H1 (1 Plate) | K049-H5 (5 Plate)

Chemiluminescent ELISA: K049-C1 (1 Plate) | K049-C5 (5 Plate)

FEATURES

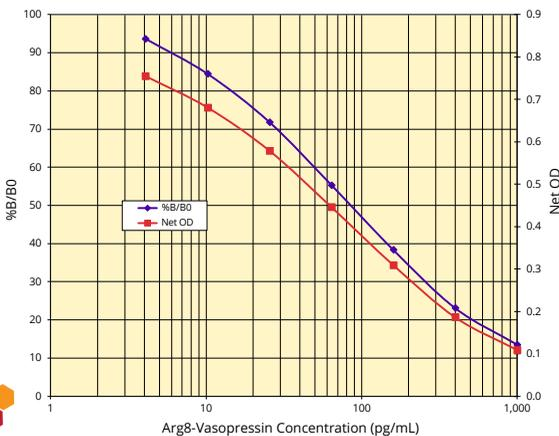
- ▶ Use: Measure AVP in Mammals, Arg-Vasotocin in Birds, Reptiles
- ▶ Sample: Extracted Serum, Plasma and Buffers, Tissue Culture Media
- ▶ Sensitivity: ELISA: 3.7 pg/mL
Chemiluminescent: < 0.9 pg/mL
- ▶ Simple: Extraction Reagent included, no C18 SPE Columns needed
- ▶ Samples/Kit: ELISA: 39 or 231 in Duplicate
Chemiluminescent: 38 or 230 in Duplicate
- ▶ Readout: ELISA: Colorimetric, 450 nm
Chemiluminescent: Glow Luminescent



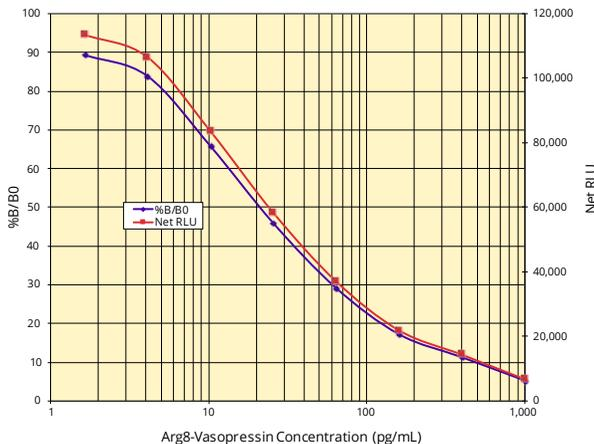
SCIENTIFIC RELEVANCE

The neurohypophysial hormone arginine vasopressin (AVP) is involved in a wide range of physiological regulatory processes, including renal water reabsorption, cardiovascular homeostasis, hormone secretion from the anterior pituitary, and modulation of social behavior and emotional status. AVP and the structurally related posterior pituitary hormone, oxytocin (OT), are synthesized in the paraventricular nucleus and the supraoptic nucleus of the hypothalamus. AVP is a 9 amino acid peptide with a 6-member disulfide ring. It is structurally related to OT, differing by 2 amino acids.

K049-H1/H5



K049-C1/C5



Ceruloplasmin (Cp) Colorimetric Activity Kit

K035-H1 (2 Plate)

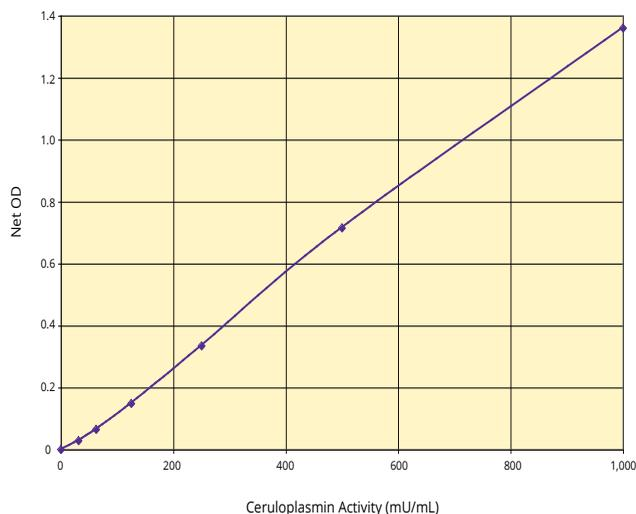
FEATURES

- ▶ 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.



Creatinine Urinary Detection Kits/Solution

Kits: K002-H1 (2 Plate) | K002-H5 (10 Plate)

Solution: X116-100ML (10 mg/dL)

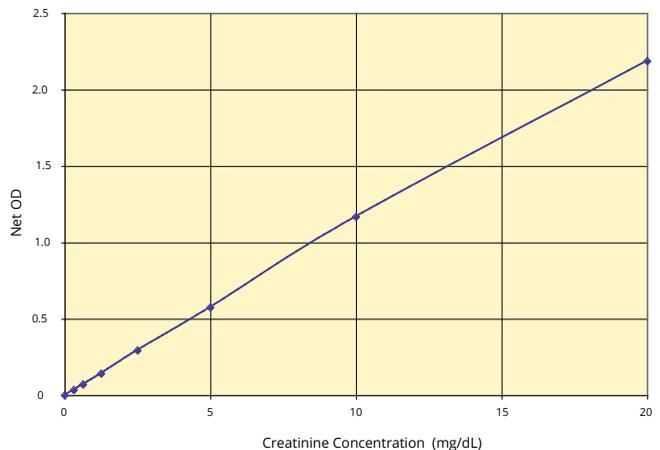
FEATURES

- ▶ Use Urine Volume Marker
- ▶ Sample Urine
- ▶ Calibrated NIST Standard Reference #914a
- ▶ Species Species Independent
- ▶ Time to Answer 30 Minutes
- ▶ Format 96-Well
- ▶ Samples/Kit 88 or 472 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Creatinine (2-amino-1-methyl-5H-imidazol-4-one) is a metabolite of phosphocreatine (p-creatinine), a molecule used as a store for high-energy phosphate that can be utilized by tissues for the production of ATP. Creatine either comes from the diet or is synthesized from the amino acids arginine, glycine, and methionine. This occurs in the kidneys and liver, although other organ systems may be involved and species-specific differences may exist. Creatine and p-creatinine are converted non-enzymatically to the metabolite creatinine, which diffuses into the blood and is excreted by the kidneys. Creatinine forms spontaneously from p-creatinine. Under normal conditions, its formation occurs at a rate that is relatively constant and as intra-individual variation is <15% from day to day, creatinine is a useful tool for normalizing the levels of other molecules found in urine. Additionally, altered creatinine levels may be associated with conditions that result in decreased renal blood flow such as diabetes and cardiovascular disease.



Dehydroepiandrosterone Sulfate (DHEA-S) ELISA Kits

K054-H1 (1 Plate) | K054-H5 (5 Plate)

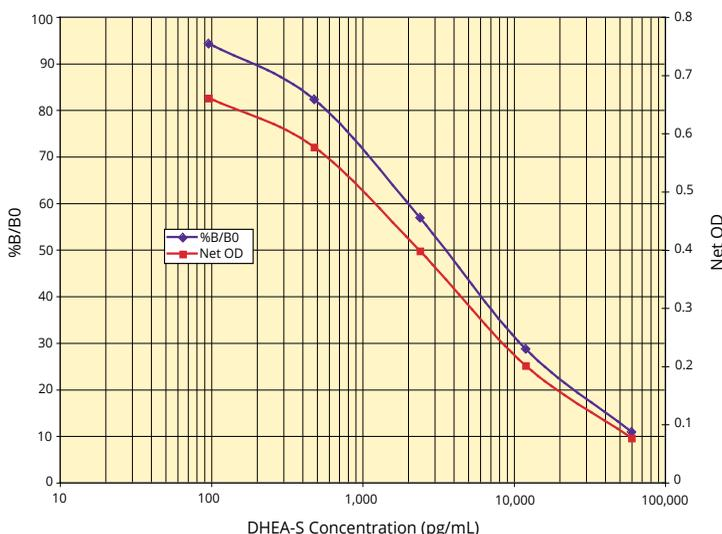
FEATURES

- ▶ Use: Determination of Estrogen Deficiency
- ▶ Sample: Serum, Plasma, Saliva, Urine, Media, Fecal Extracts
- ▶ Validation: Human, Monkey, Felids, Ungulates
- ▶ Time to Answer: 2.5 Hours
- ▶ Format: 96-Well, Break-Apart Strip
- ▶ Species: Species Independent
- ▶ Samples/Kit: 41 or 233 in Duplicate
- ▶ Stability: Liquid 4°C Stable Reagents
- ▶ Readout: Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Dehydroepiandrosterone sulfate (DHEA-S) is the major C19 steroid secreted by the adrenal cortex, and is a precursor to testosterone and estrogen biosynthesis. It is produced by the addition of a sulfate group to dehydroepiandrosterone (DHEA) catalyzed by the sulfotransferase enzymes, SULT1A1 and SULT1E1, which also produce estrone sulfate from estrone. Due to the presence of a 17-ketone group rather than a hydroxyl group, DHEA-S has relatively low androgenic activity. The bioactivity of DHEA-S may be high due to its serum concentrations being 100-1,000-fold higher than testosterone or DHEA in addition to its weak affinity for sex-hormone binding globulin.



Epiandrosterone ELISA Kits

K063-H1 (1 Plate) | K063-H5 (5 Plate)

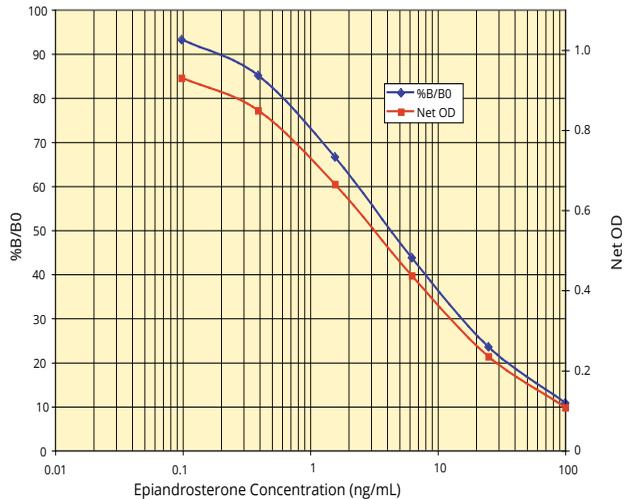
FEATURES

- ▶ Use Measure Epiandrosterone in a Variety of Matrices
- ▶ Sample Fecal Extracts, Urine, Saliva, Extracted Serum/Plasma
- ▶ Sensitivity 0.120 ng/mL
- ▶ Time to Answer 2.5 Hours
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Epiandrosterone is a naturally occurring metabolite of dehydroepiandrosterone (DHEA) found in most mammals and produced via the action of the 5 α -reductase enzyme. It is a weak androgen formed primarily in peripheral tissues, released into circulation and ultimately excreted in urine. Epiandrosterone has been shown to inhibit the pentose phosphate pathway (PPP), decreasing intracellular NADPH levels. It also attenuates NO-evoked relaxation of the pulmonary artery. It has been linked to gonadal activity and sexual behavior in males. Epiandrosterone is of interest for cell metabolism, cardiac, and prostate cancer research.

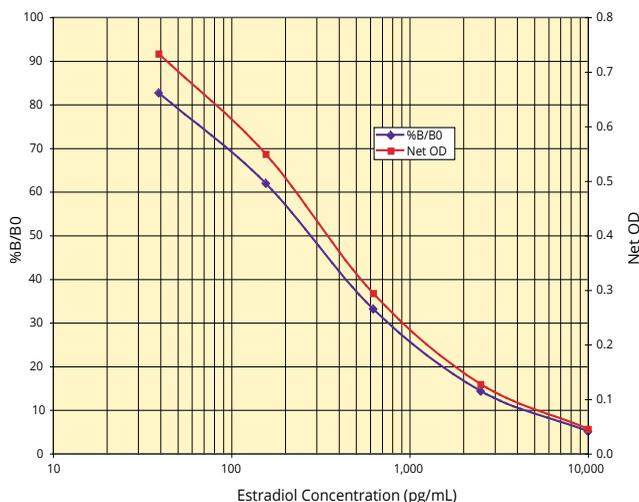


FEATURES

- ▶ Use Noninvasive Estrogen Assessment
- ▶ Sample Urine, Media, and Fecal Extracts
- ▶ Validation Mice, Rats, Humans, Monkeys, Birds, Felids, Ungulates
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 41 or 233 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm

SCIENTIFIC RELEVANCE

Estradiol (E2, 17 β -estradiol, or oestradiol) is the predominant sex hormone present in females. It is also present in males, being produced as an active metabolic product of testosterone. It represents the major estrogen in humans. Estradiol has not only a critical impact on reproductive and sexual functioning, but also affects other organs. Serum estradiol measurements in women reflect primarily the activity of the ovaries. As such, they are useful in the detection of baseline estrogen in women with amenorrhea or menstrual dysfunction and to detect the state of hypoestrogenism and menopause. Furthermore, estrogen monitoring during fertility therapy assesses follicular growth. Estrogen-producing tumors and in precocious puberty samples will demonstrate persistent high levels of estradiol and other estrogens.



Serum Estradiol ELISA Kits

KB30-H1 (1 Plate) | KB30-H5 (5 Plate)

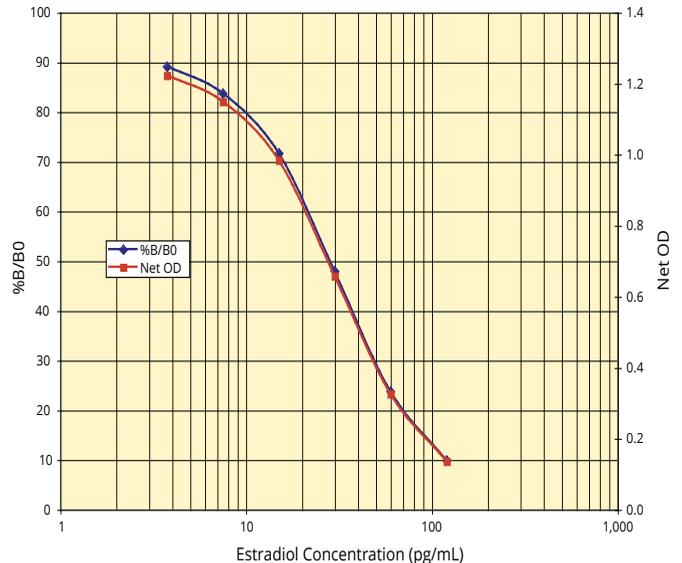
FEATURES

- ▶ Use Estrogen Assessment
- ▶ Sample Serum and Plasma
- ▶ Validation Mice, Rats, Humans, Monkeys, Birds, Felids, Ungulates
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 for in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Estradiol (E2, 17 β -estradiol, or oestradiol) is the predominant sex hormone present in females. It is also present in males, being produced as an active metabolic product of testosterone. It represents the major estrogen in humans. Estradiol has not only a critical impact on reproductive and sexual functioning, but also affects other organs. Serum estradiol measurements in women reflect primarily the activity of the ovaries. As such, they are useful in the detection of baseline estrogen in women with amenorrhea or menstrual dysfunction and to detect the state of hypoestrogenism and menopause. Furthermore, estrogen monitoring during fertility therapy assesses follicular growth. Estrogen-producing tumors and in precocious puberty samples will demonstrate persistent high levels of estradiol and other estrogens.



FEATURES

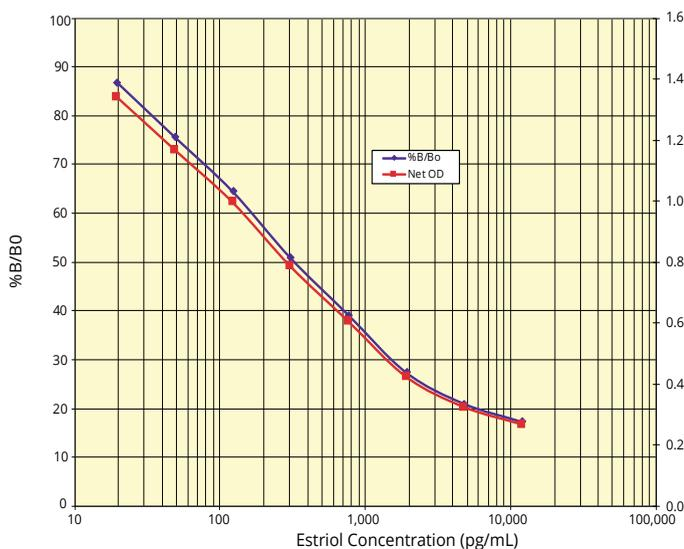
- ▶ Use Estrogen Assessment
- ▶ Sample Urine, Saliva, and Fecal Extracts
- ▶ Time to Answer 2.5 Hours or Overnight
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 38 or 230 in Duplicate
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Estriol is one of the three major endogenous estrogens, along with estradiol and estrone. It is a weak estrogen derived from hydroxylation of estradiol and estrone in the liver. Normal levels in women who are not pregnant are typically nearly undetectable. However, estriol is produced in large amounts by the placenta and rising maternal levels can be detected from the very early weeks of pregnancy through until delivery.

Estriol can be monitored as an indicator of fetal health and well-being during pregnancy. It is routinely measured as part of both the triple test and the quadruple test during pregnancy outreach and screening. Abnormally low levels in pregnant females can suggest chromosomal or congenital anomalies in the fetus. In some parts of the world exogenous estriol is used for the treatment of menopausal symptoms. Estriol has also been investigated as a protective neurosteroid with potential roles in immune diseases and bone and lipid metabolism.



Estrone ELISA Kits

K031-H1 (1 Plate) | K031-H5 (5 Plate)

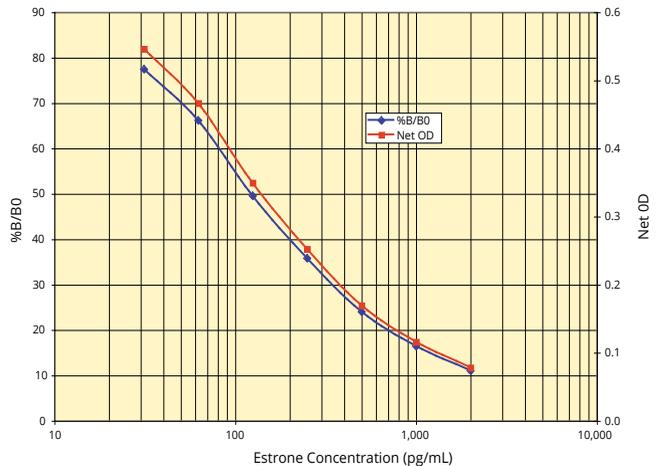
FEATURES

- ▶ Use Determination of Inborn Errors of Sex Steroid Metabolism
- ▶ Sample Urine, Fecal Extracts, and Media
- ▶ Validation Multiple Species
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 39 or 231 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Estrone, also known as E1 or oestrone (3-hydroxy-1,3,5(10)-estratrien-17-one) is a C-18 steroid hormone and is one of the three naturally occurring estrogens, the others being estradiol and estriol. Estrone is produced primarily from androstenedione, originating from the gonads or the adrenal cortex and from estradiol by 17-hydroxysteroid dehydrogenase enzyme systems. Estrone concentrations in premenopausal mammals fluctuate according to the menstrual cycle. In premenopausal women, more than 50% of the estrone is secreted by the ovaries. Interconversion of estrone and estradiol also occurs in peripheral tissue. In humans, during the follicular phase of the menstrual cycle estrone levels increase slightly.



Estrone-3-Glucuronide (E1G) ELISA Kits

K036-H1 (1 Plate) | K036-H5 (5 Plate)

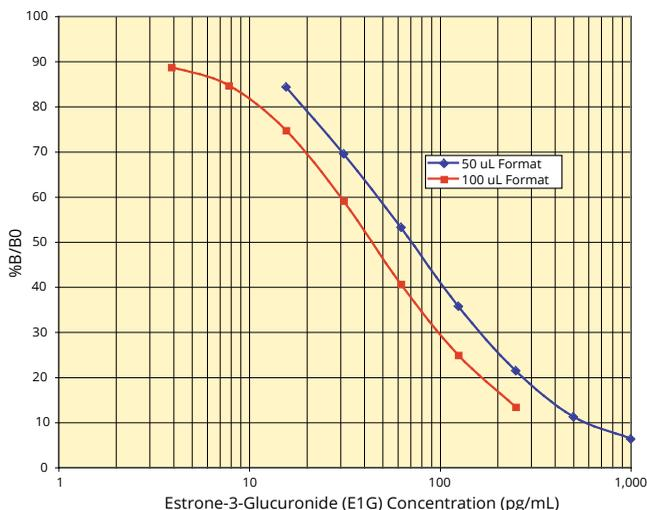
FEATURES

- ▶ Use Estrogen Assessment
- ▶ Sample Urine, Media, and Fecal Extracts
- ▶ Validation Multiple Species
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 39 or 231 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Estrone-3-glucuronide (1,3,5(10)-estratrien-3-ol-17-one glucosiduronate, E1G) is the principle secreted form of circulating estradiol in mammals. Ovulation is the critical event of each menstrual cycle that occurs during the reproductive life of healthy females and the ovum can only be fertilized during the short period of time in which it is viable. The three phases of the menstrual cycle are: (i) an initial phase when there is only a low risk that would enable viable spermatozoa to survive and reach the ovum, (ii) a phase when the chance of fertilization is at a maximum, the fertile period, and (iii) a time of absolute infertility when the ovum is no longer viable. Clinical studies have indicated the utility of measuring estrone-3-glucuronide (E1G) and pregnanediol-3-glucuronide (PDG) in samples of urine or fecal extracts to monitor ovarian function in females.



Estrone-3-Sulfate (E1S) ELISA Kits

K038-H1 (1 Plate) | K038-H5 (5 Plate)

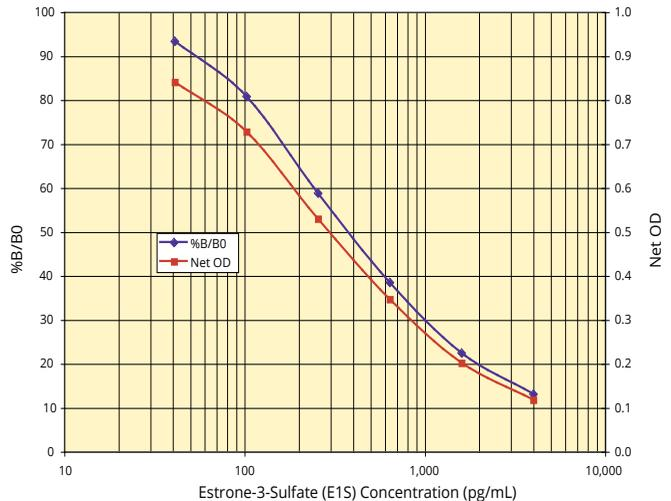
FEATURES

- ▶ Use Breast Cancer and Cryptorchidism
- ▶ Sample Serum, Plasma, Urine, Fecal Extracts and Media
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Estrone-3-sulfate (E1S) is synthesized in the fetal or cotyledonary portion of the placenta. Estrone sulfate, which is present in plasma at a higher concentration than either unconjugated estrone or estradiol in nonpregnant women and normal men, appears to originate almost entirely from a conjugation of estrone and converted estradiol in non-glandular tissues. Estrone sulfate is quantitatively the most important circulating estrogen. Breast tumors contain sulfatase activity and can convert estrone sulfate into estradiol. Cryptorchidism, where one or both testicles fail to descend is considered to be a prevalent defect in horses. Bilaterally cryptorchid stallions do not produce viable spermatozoa but often exhibit normal secondary sexual characteristics. Several investigators have suggested measuring testosterone and estrone sulfate serum levels as reliable diagnostic aids for the condition.



Levonorgestrel (LNG) ELISA Kits

K058-H1 (1 Plate) | K058-H5 (5 Plate)

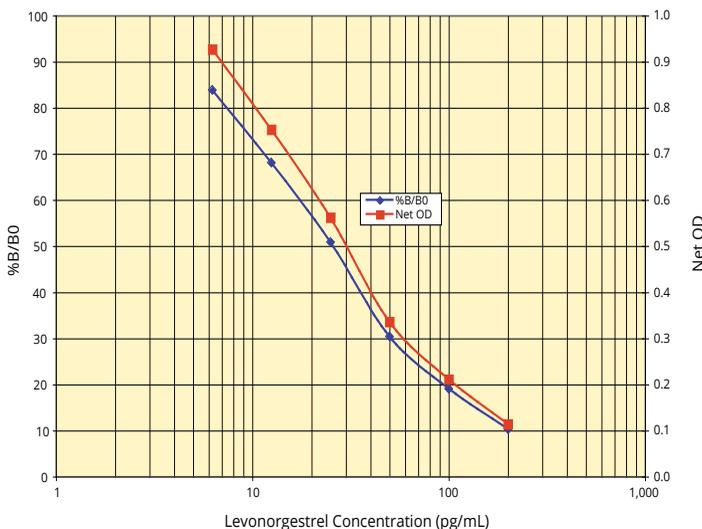
FEATURES

- ▶ Use Measure LNG in a Variety of Matrices
- ▶ Sample Saliva, Urine, Water, Milk, TCM, and Extracted Serum, Plasma and Fecal Material
- ▶ Time to Answer 1.5 Hours
- ▶ Sensitivity 2.20 pg/mL
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Levonorgestrel (LNG) is a synthetic steroid commonly used for contraception, treatment of dysmenorrhea, and for endometrial protection during estrogen replacement therapy in postmenopausal women. LNG has also been shown to be an effective treatment modality for a variety of gynecologic conditions including: heavy menstrual bleeding, pelvic pain, endometrial hyperplasia and early stage endometrial cancer. LNG is safe to use while breastfeeding and works by decreasing ovulation in addition to changing the mucus in the cervix to prevent the passage of sperm and altering the uterine lining. Quantitative measurement of LNG in biological samples can be useful for determining if target therapeutic concentrations are being met and maintained. LNG can also be measured in environmental samples using this kit. Environmental LNG can have toxic effects in aquatic ecosystems.



Oxytocin ELISA & Chemiluminescent ELISA Kits/Solutions

ELISA: K048-H1 (1 Strip-Well Plate) | K048-H5 (5 Strip-Well Plate)

ELISA: K048-H1W (1 Whole Plate) | K048-HSW (5 Solid Plate)

Chemiluminescent ELISA: K048-C1 (1 Plate) | K048-C5 (5 Plate)

Solutions: Isotocin X128-625UL | Mesotocin X127-625UL

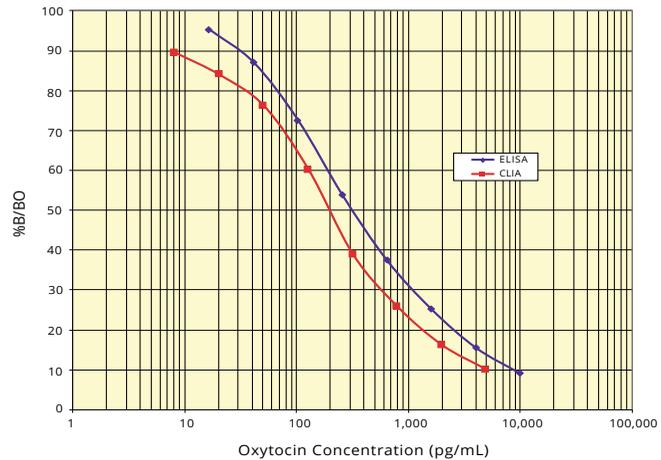


FEATURES

- ▶ Use Reproductive Assessment
- ▶ Sample Clarified Milk and Extracted Serum, Plasma, Saliva
- ▶ Time to Answer Overnight
- ▶ Sensitivity ELISA: 17.0 pg/mL
Chemiluminescent: 6.33 pg/mL
- ▶ Species Mammals, Birds, Fish
- ▶ Cross Reactivity High Reactivity to Mesotocin (88.4%) and Isotocin (94.3%)
- ▶ Samples/Kit 38 or 230 in Duplicate
- ▶ Readout ELISA: Colorimetric, 450 nm
Chemiluminescent: Glow Luminescence

SCIENTIFIC RELEVANCE

The neuropeptides oxytocin and vasopressin were isolated and synthesized by Vincent du Vigneaud, work for which he received the Nobel Prize in Chemistry in 1955. Oxytocin is a neurohypophysial peptide produced in the hypothalamus. The molecule consists of nine amino acids linked with a [1-6] disulfide bond and a semi-flexible carboxyamidated tail. Highly conserved across species boundaries, oxytocin-like neurohypophysial peptides are substituted primarily at residues 4 and/or 8. In the oxytocin-like peptide, mesotocin; a common peptide found in some fishes, reptiles, amphibians, marsupials and non-mammalian tetrapods, the leucine at residue 8 is substituted for isoleucine. Oxytocin binds to specific cell surface receptors, which in turn initiate a secondary intracellular response cascade via a phosphoinositide signaling pathway.



PGFM (13,14-Dihydro-15-keto-Prostaglandin F2) ELISA Kits

K022-H1 (1 Plate) | K022-H5 (5 Plate)

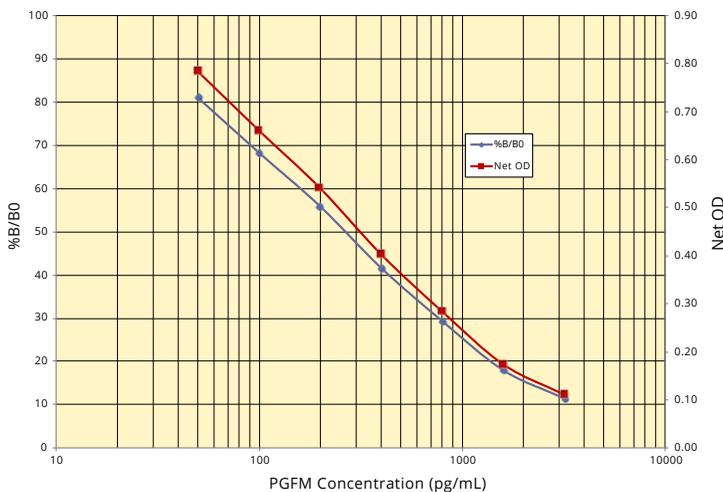
FEATURES

- ▶ Use: Pregnancy Marker
- ▶ Sample: Serum, Plasma, Urine, Fecal Extracts and Media
- ▶ Validation: Felids and Pandas
- ▶ Time to Answer: 90 Minutes
- ▶ Format: 96-Well, Break-Apart Strip
- ▶ Species: Species Independent
- ▶ Samples/Kit: 39 or 231 in Duplicate
- ▶ Stability: Liquid 4°C Stable Reagents
- ▶ Readout: Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Uterine and placental Prostaglandin F₂ (PGF₂) is involved in the regulation of reproductive and pregnancy-related processes such as embryonic development, initiation of parturition, and resumption of ovarian activity. In domestic ruminants, uterine tissue is a primary source of PGF₂ and secretion of uterine PGF₂ is a key regulator for the cyclical regression of the corpus luteum. Prostaglandin F₂ is metabolized to PGFM (13,14-dihydro-15-keto-PGF₂) during passage through the lungs. PGFM has a longer half-life in peripheral circulation than PGF₂ and has been applied as a useful analytical marker of PGF₂. PGFM is a useful non-invasive marker of pregnancy when measured in both urine and fecal samples. It has been shown to be a precise, practical method for this application in these matrices. Fecal PGFM analyses may allow pregnancy diagnosis in captive and free-ranging felids as well as pandas and other species.



Pregnanediol-3-Glucuronide (PDG) ELISA Kits

K037-H1 (1 Plate) | K037-H5 (5 Plate)

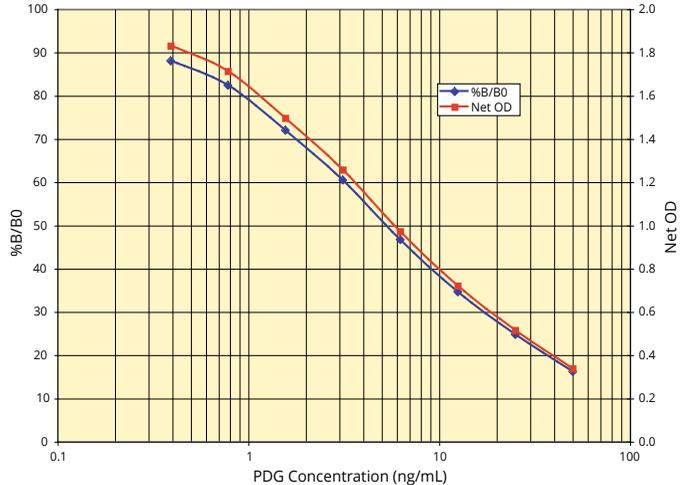
FEATURES

- ▶ Use Pregnancy Assessment
- ▶ Sample Urine, Media and Extracted Serum, Plasma, Fecal
- ▶ Validation Multiple Species
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Species Species Independent
- ▶ Samples/Kit 41 or 233 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm

**MULTI
SPECIES**

SCIENTIFIC RELEVANCE

Pregnanediol-3-Glucuronide also known as PDG 5 β -pregnane-3 α ,20 α -diol 3 α -glucuronide; is the major metabolite of progesterone. Progesterone is a hormone involved in the female menstrual cycle, gestation and embryogenesis of humans and other species. Progesterone belongs to a class of hormones called progestogens and is the major naturally occurring human progestogen. Progesterone is an essential regulator of human female reproductive function in the uterus, ovary, mammary gland and brain, and play important roles in the cardiovascular, skeletal, and central nervous systems. Progesterone also has neurotrophic roles in the peripheral nervous system as it activates the growth and maturation of axons and stimulates the repair and replacement of myelin sheaths in regenerating nerve fibres.



Progesterone ELISA Kits

K025-H1 (1 Strip-Well Plate) | K025-H5 (5 Strip-Well Plate)
 K025-H1W (1 Whole Plate) | K025-H5W (5 Whole Plate)

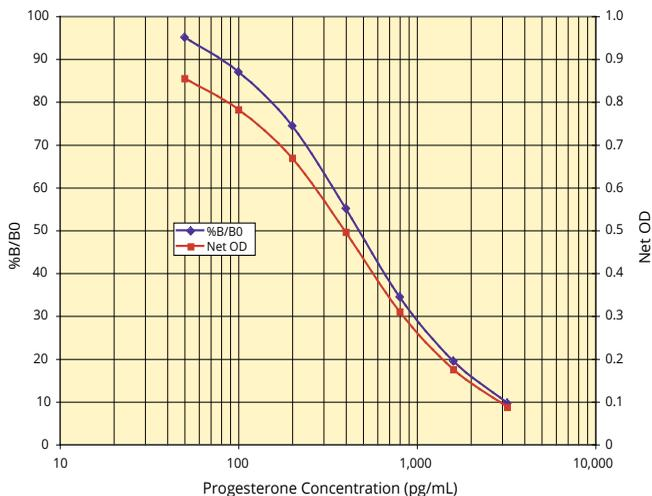


FEATURES

- ▶ Use: Determination of Reproduction and Sex Steroid Metabolism
- ▶ Sample: Urine, Fecal Extracts, Media, Serum and Plasma
- ▶ Validation: Multiple Species
- ▶ Time to Answer: 2.5 Hours
- ▶ Sensitivity: 47.9 pg/mL
- ▶ Species: Species Independent
- ▶ Samples/Kit: 39 or 231 in Duplicate
- ▶ Stability: Liquid 4°C Stable Reagents
- ▶ Readout: Colorimetric, 450 nm

SCIENTIFIC RELEVANCE

Progesterone, also known as P4 (Pregn-4-ene-3,20-dione), is a C21 steroid hormone involved in the female menstrual cycle, gestation and embryogenesis of humans and other species. Progesterone belongs to a class of hormones called progestogens and is the major naturally occurring progestogen. Progesterone is an essential regulator of human female reproductive function in the uterus, ovary, mammary gland and brain, and plays important roles in non-reproductive tissues such as the cardiovascular, skeletal, and central nervous systems. Progesterone action is conveyed by two isoforms of the nuclear progesterone receptor (PR), PR-A and PR-B. PR-A and -B are expressed in a variety of normal breast tissue from humans, rats and mice and is also expressed in breast cancer cells. Progesterone also has neurotrophic roles in the peripheral nervous system as it activates the growth and maturation of axons and stimulates the repair and replacement of myelin sheaths in regenerating nerve fibres.



Progesterone Metabolites ELISA Kits

K068-H1 (1 Plate) | K068-H5 (5 Plate)

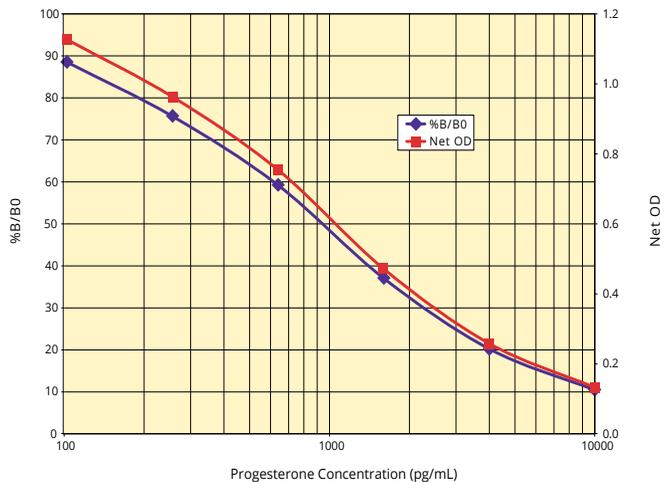
FEATURES

- ▶ Use Measure general progesterone metabolites and generate reproductive profiles
- ▶ Sample Fecal Extracts, Urine
- ▶ Time to Answer 90 Minutes
- ▶ Sensitivity 51.2 pg/mL
- ▶ Species Species Independent
- ▶ Samples/Kit 40 or 232 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Progesterone (P4, pregn-4-ene-3,20-dione) belongs to a class of hormones called progestogens, and is an essential regulator of female reproductive function in the uterus, ovary, mammary gland and brain. Progestogens are the primary hormones involved in the female menstrual cycle, gestation, and embryogenesis of humans and most other species. Progesterone can be metabolized and excreted as a variety of general progesterone by-products. Common metabolites include 5-reduced progesterone (pregnane), pregnanolones and hydroxyprogesterones. Measurement of progesterone metabolites provides vital data about reproductive status and is essential for studying reproductive and survival strategies of wildlife and endangered species. Likewise, assays measuring progesterone metabolites provide a clearer picture of hormonal regulation and reproduction.



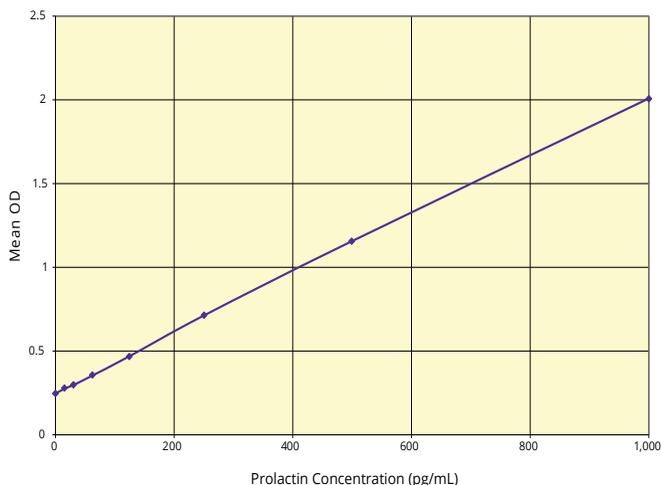
FEATURES

- ▶ Use Reproduction
- ▶ Sample Serum and Plasma
- ▶ Validation Humans and Elephants
- ▶ Time to Answer 2.5 Hours
- ▶ Format 96-Well, Break-Apart Strip
- ▶ Samples/Kit 40 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Prolactin (PRL) is a polypeptide hormone that is synthesized and secreted from specialized cells of the anterior pituitary gland. The hormone was given its name based on the fact that an extract of bovine pituitary gland would cause growth of the crop sac and stimulate the production of milk in pigeons or promote lactation in rabbit. However, it is now appreciated that prolactin has over 300 separate biological activities. Prolactin has multiple roles in reproduction other than lactation, and it also plays multiple homeostatic roles in the organism. Furthermore, the synthesis and secretion of prolactin is not restricted to the anterior pituitary gland, but multiple other organs and tissues in the body have this capability.



Testosterone ELISA Kits

Strip Well: K032-H1 (1 Plate) | K032-H5 (5 Plate)

Whole Plate: K032-H1W (1 Plate) | K032-H5W (5 Plate)

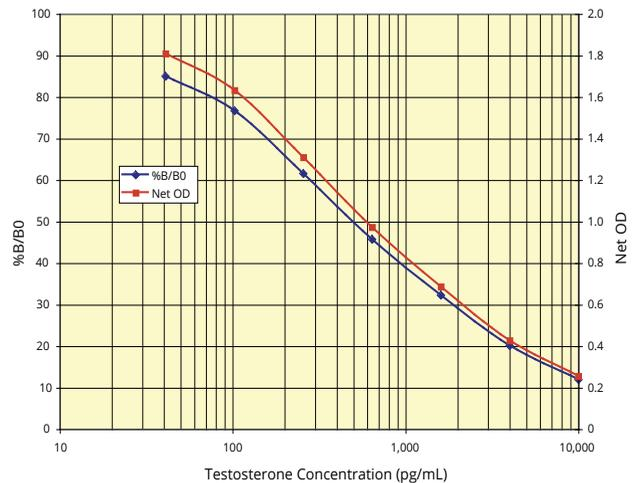
FEATURES

- ▶ Use Non-invasive measurement of Testosterone and Dihydrotestosterone
- ▶ Sample Dried Fecal Extracts, Urine, TCM, Extracted Serum and Plasma
- ▶ Range 40.96- 10,000 pg/mL
- ▶ Sensitivity 9.92 pg/mL
- ▶ Time to Answer 2.5 Hours
- ▶ Species Species Independent
- ▶ Samples/Kit 39 or 231 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Testosterone (4-Androsten-17 β -ol-3-one) is an anabolic steroid hormone from the androgen group. It is found in mammals, reptiles, birds, and other vertebrates. In mammals, testosterone is primarily secreted in the testes of males and the ovaries of females, although small amounts are also secreted by the adrenal glands. It is the principal male sex hormone and plays key roles in the development of reproductive tissues such as the testes and prostate, and in promoting secondary sexual characteristics such as increased muscle, bone mass, and body hair. In addition, testosterone is essential for health and well-being as well as the prevention of osteoporosis. Testosterone plays a significant role in glucose homeostasis and lipid metabolism. Cross-sectional epidemiological studies have reported a direct correlation between plasma testosterone and insulin sensitivity. Low testosterone levels are associated with an increased risk of type 2 diabetes, dramatically illustrated by androgen deprivation in men with prostate carcinoma.



Testosterone ELISA Kits - Improved Sensitivity

Strip-Well: K080-H1 (1 Plate) | K080-H5 (5 Plate)

Whole Plate: K080-H1W (1 Plate) | K080-H5W (5 Plate)

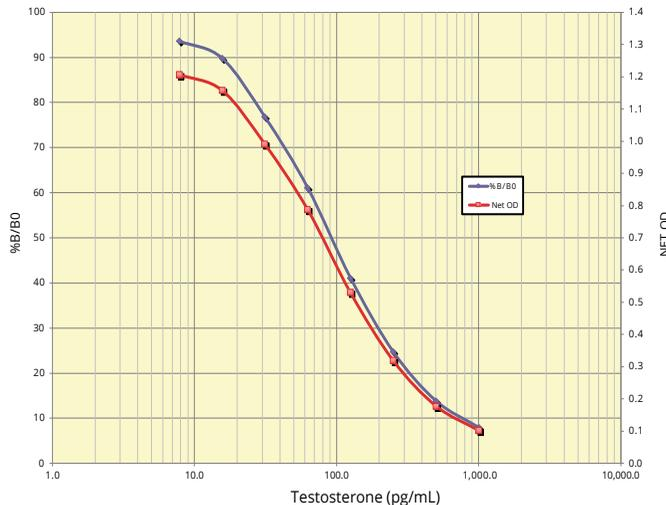
FEATURES

- ▶ Use No Extraction Needed for Serum and Plasma
- ▶ Sample Serum, Plasma, Urine and Dried Fecal Extracts
- ▶ Range 7.813 -1,000 pg/mL
- ▶ Sensitivity 2.97 pg/mL
- ▶ Time to Answer 2.5 Hours
- ▶ Species Species Independent
- ▶ Samples/Kit 38 or 230 in Duplicate
- ▶ Stability Liquid 4°C Stable Reagents
- ▶ Readout Colorimetric, 450 nm



SCIENTIFIC RELEVANCE

Testosterone (4-Androsten-17β-ol-3-one) is an anabolic steroid hormone from the androgen group. It is found in mammals, reptiles, birds, and other vertebrates. In mammals, testosterone is primarily secreted in the testes of males and the ovaries of females, although small amounts are also secreted by the adrenal glands. It is the principal male sex hormone and plays key roles in the development of reproductive tissues such as the testes and prostate, and in promoting secondary sexual characteristics such as increased muscle, bone mass, and body hair. In addition, testosterone is essential for health and well-being as well as the prevention of osteoporosis. Testosterone plays a significant role in glucose homeostasis and lipid metabolism. Cross-sectional epidemiological studies have reported a direct correlation between plasma testosterone and insulin sensitivity. Low testosterone levels are associated with an increased risk of type 2 diabetes, dramatically illustrated by androgen deprivation in men with prostate carcinoma.





STRESS

REPRODUCTIVE

OXIDATIVE STRESS



METABOLISM

VOLUME & NORMALIZATION



KIDNEY INJURY



INFLAMMATION



CELL SIGNALING



ARBOR ASSAYS™