



Wildlife Endocrinology

Steroid Hormone ELISA Kits

Enzo's Trusted Steroid Hormone ELISA Kits Deliver Sensitivity and Reproducibility

With thousands of species each year listed as endangered or threatened, the work done by biologists around the world to understand and preserve animal health is never ending. Reproduction studies and stress are two of the most common concerns today. Hormone analysis continues to be widely accepted for monitoring stress response and evaluating reproductive health on wildlife. Enzo Life Sciences provides a broad range of immunoassay kits for monitoring stress, reproductive physiology, health and welfare, and survival of wildlife animals.

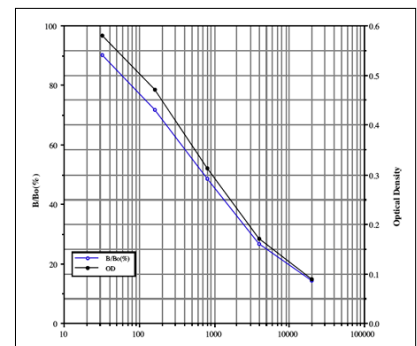
STRESS MONITORING

Corticosterone ELISA Kit (ADI-900-097)

- Sensitive measurement of corticosterone, detecting as little as 27 pg/mL
- High-throughput with results for up to 39 samples in duplicate in just 3 hours
- Built-in flexibility to measure corticosterone in any species
- Trusted and routinely used to monitor stress levels in zoo animals

PRODUCT	PRODUCT #	VALIDATED SAMPLES	SPECIES
Aldosterone ELISA Kit	ADI-900-173	P, S, U	Species Independent
Angiotensin II ELISA Kit	ADI-900-204	P, S	Species Independent
Corticosterone ELISA Kit	ADI-900-097	CS, P, S, Sa, and F	Species Independent
Cortisol ELISA KIT	ADI-900-071	CS, P, S, Sa, U	Species Independent
Direct 8-iso-PGF ₂ α ELISA Kit	ADI-900-091	P, S, TL	Species Independent

Highly Sensitive Measurement Of Corticosterone



Typical standard curve of Corticosterone ELISA Kit (ADI-900-097)

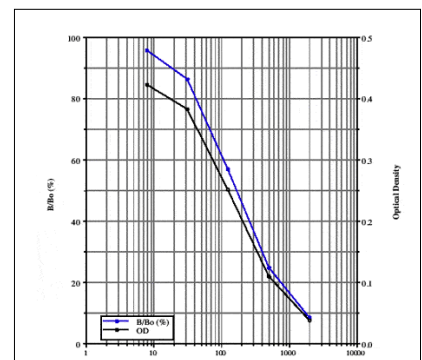
REPRODUCTIVE STUDIES

Testosterone ELISA Kit (ADI-900-065)

- Highly sensitive measurement of testosterone, detecting as little as 5.67 pg/mL
- Steroid displacement reagent provided to assay free or total testosterone
- High-throughput format with results in 3 hours for up to 39 samples in duplicate
- Excellent lot-after-lot consistency ideal for longitudinal studies

PRODUCT	PRODUCT #	VALIDATED SAMPLES	SPECIES
17β-Estradiol ELISA Kit	ADI-900-008	CS, Sa	Species Independent
17β-Estradiol High Sensitivity ELISA Kit	ADI-900-174	P, S	Species Independent
Arg ⁸ -Vasopressin ELISA Kit	ADI-900-017A	CS, P, S	Species Independent
DHEA ELISA Kit	ADI-900-093	CS, P, S, Sa, U	Species Independent
Direct 8-iso-PGF ₂ α ELISA Kit	ADI-900-091	P, S, TL	Species Independent
Oxytocin ELISA Kit	ADI-900-153A	CS, S, Sa, U, CSF	Species Independent
Progesterone ELISA Kit	ADI-900-011	CS, S, Sa	Species Independent
Testosterone ELISA Kit	ADI-900-065	CS, P, S, Sa	Species Independent
Testosterone High Sensitivity ELISA Kit	ADI-900-176	P, S, U	Species Independent

Highly Sensitive Measurement Of Testosterone



Typical standard curve of Testosterone ELISA Kit (ADI-900-065)

KEY: CS = Culture supernatant; CSF = Cerebral Spinal Fluid; F = Feces; P = Plasma; S = Serum; Sa = Saliva; TL = Tissue Lysate; U = Urine

Advance Your Research:

www.enzolifesciences.com/wildlife-endocrinology

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com

Rev. 082420