## Oxytocin ELISA kit

The most sensitive, specific and flexible alternative to more costly and time-consuming methods for oxytocin detection.

The Oxytocin EIA kit is a colorimetric competitive enzyme immunoassay kit with results overnight + 1 hour. Absorbance is read at 405 nm. Very low reactivity with vasopressin provides confidence in assay results. Sensitive quantification of 15pg/mL oxytocin. Ready-to-use liquid color-coded reagents reduce errors in the lab.

Oxytocin is a neurohypophysial peptide which is produced in the paraventricular nuclei of the hypothalamus and stored in the posterior pituitary. The molecule consists of nine amino acids linked with a [1-6] disulfide bond and a semi-flexible carboxyamidated tail. A hormone once thought to be limited to female smooth muscle reproductive physiology, more current findings have determined that oxytocin also functions as a neurotransmitter, may be involved in neuropsychiatric disorders, social/sexual behavior, and is important in male reproductive physiology. Oxytocin and the related neurohypophysial peptide, Arg-Vasopressin, maintain renal water and sodium balance.

Citations: 185

View Online »

**Ordering Information** 

Order Online »

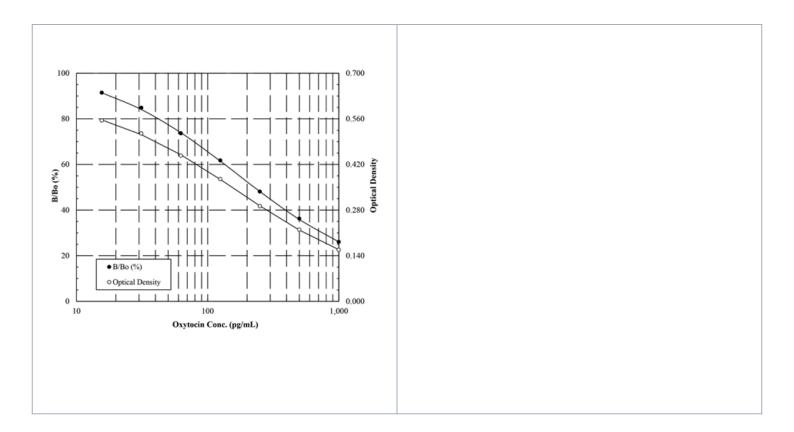
ADI-900-153A-0001

96 wells

Manuals, SDS & CofA

View Online »

- Highly specific ELISA which does not detect vasopressin, providing confidence in assay results
- Ultra-sensitive measurement, detecting as little as 15pg/ml oxytocin
- Cost effective solution, faster and less costly than LC/MS
- Easy-to-use liquid color-coded reagents and pre-coated plates minimizes errors
- Widely published in peer reviewed literature



## **Handling & Storage**

**Use/Stability** Store all components at +4°C, except standard and conjugate at -20°C.

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application** Colorimetric detection, ELISA

**Application Notes**For the quantitative determination of Oxytocin in culture

supernatants, milk, plasma, and serum from any species. Validated sample types include cerebral spinal fluid, saliva,

tissue and urine.

Assay Time Overnight + 1 hour

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

**Contents** GxR IgG Microtiter plate, Conjugate, Antibody, Assay

buffer, Wash buffer concentrate, Standard, pNpp

Substrate, Stop solution

Crossreactivity Mesotocin (7%), Oxytocin (100%), Arg8-Vasotocin (7.5%),

and <0.02% for other related molecules.

**Sensitivity** 15pg/ml (range 15.6 – 1,000 pg/ml)

Species Reactivity Species independent

Technical Info / Product Notes Cited samples:

Hormone ELISAs: Cited Samples

Wildlife Endocrinology: Cited Sample Types

The Oxytocin ELISA kit has been extensively benchmarked for behavioral studies with plasma samples from a variety of species. **Featured in:** Frontiers in Neuroscience – Validation of a commercially available enzyme immunoassay for the determination of oxytocin in plasma samples from seven domestic animal species.

UniProt ID P01178 (human)

Wavelength 405 nm

Last modified: May 29, 2024

