

# 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 carboxyamided 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:** 186

[View Online »](#)

## Ordering Information

[Order Online »](#)

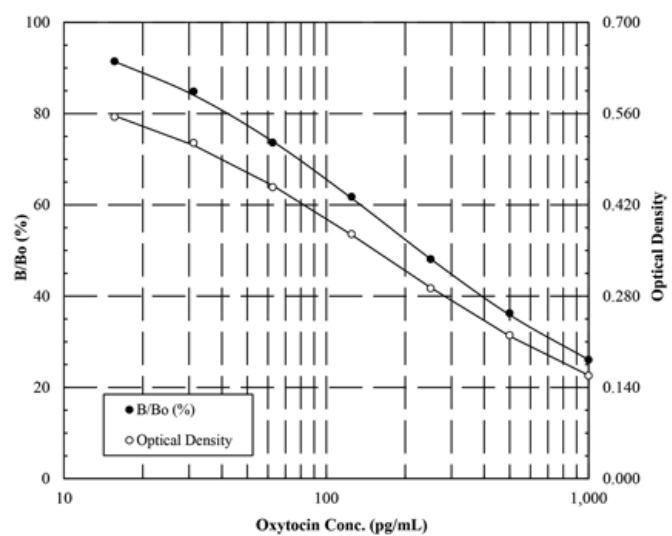
ADI-901-153A-0001

5x96 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 15 pg/mL oxytocin
- Cost effective solution, faster and less costly than LC/MS
- Easy-to-use liquid color-coded reagents and pre-coated plates minimize errors
- Widely published in peer reviewed literature
- Save with bulk kit package!



## 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 Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, 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



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)