

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: 186

[View Online »](#)

Ordering Information

[Order Online »](#)

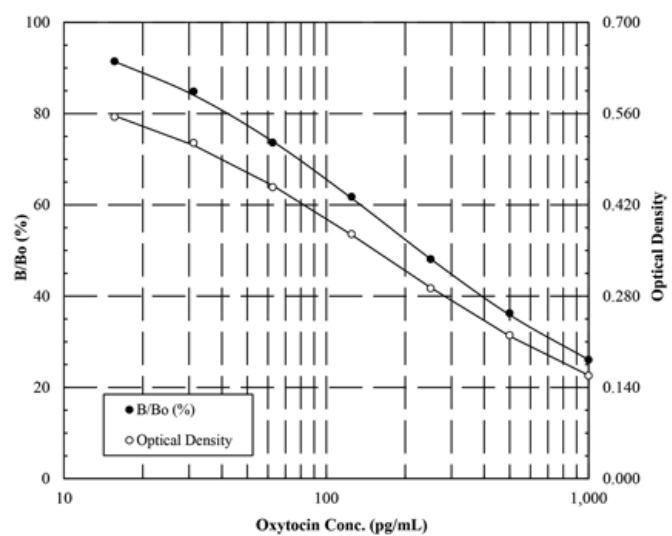
ADI-900-153A-0001

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



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)