

# Serotonin ELISA kit

Easy-to-use, well-validated serotonin ELISA kit for hormone research.

The Serotonin EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours.

Serotonin (5-hydroxytryptamine, 5-HT) is a monoamine found in the central nervous system, gastrointestinal tract, and blood with broad physiological functions in neurotransmission, gastric motility, hemostasis, and cardiovascular integrity. Seven subclasses (5-HT 1-7) of serotonin receptors have been identified, a majority of which are members of the seven transmembrane domain-containing GPCR family, and are the target of several anti-depressant drugs. The 5-HT<sub>1</sub> subclass consists of five highly homologous receptors that couple via G(i/o) alpha subunits to inhibit adenylyl cyclase activity. 5-HT<sub>1A</sub> expression in mammals is strongest in CNS limbic pathways that control emotion, including tissue of the hippocampus, dorsal raphe, and septum. Serotonin is a key biomarker used for the monitoring of overall welfare and/or the assessment of behavioral disorders in a variety of species.

Citations: 105

[View Online »](#)

## Ordering Information

[Order Online »](#)

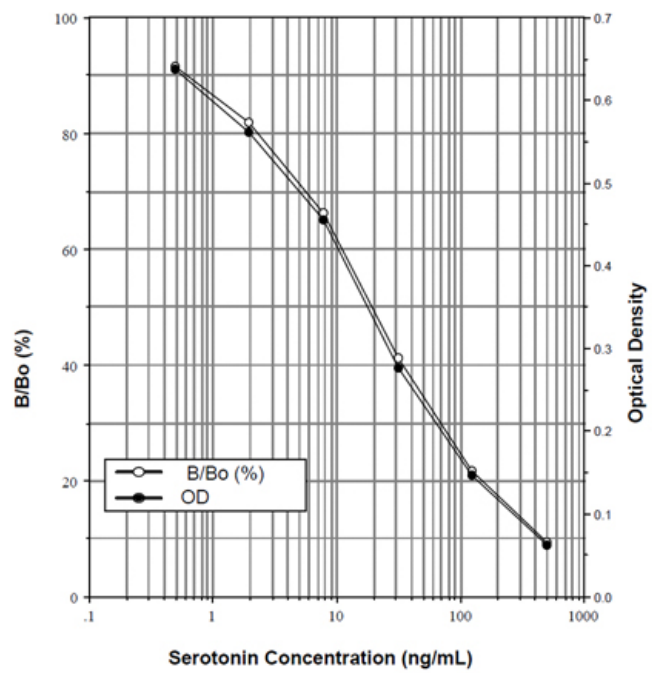
ADI-900-175

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement of serotonin, detecting as little as 0.293 ng/ml
- No acylation required for standards and samples
- Thoroughly validated in multiple complex sample matrices
- High throughput format with results in just 3 hours for up to 39 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis



## Handling & Storage

**Use/Stability** Store all components at -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5-HT, 5-hydroxytryptamine

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of Serotonin in platelets, plasma, serum, and urine from any species. Cited sample type includes culture supernatant.

**Assay Time** 3 hours

**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** N-acetyl serotonin (17%), 5-hydroxy-L-tryptophan (0.4%), Tryptamine (0.1%), 5-hydroxyindoleacetic acid (0.03%), Melatonin (0.01%), Tyramine (<0.004%), Tryptophan (<0.004%)

**Sensitivity** 0.293 ng/ml (range 0.49 – 500 ng/ml)

**Species Reactivity** Species independent

**Technical Info / Product Notes** **Application Note:**  
[Validation of a Serotonin ELISA Kit with Blood Samples from Three Domestic Animal Species.](#)

**Cited samples:**  
[Hormone ELISAs: Cited Samples](#)

**Wavelength** 405 nm



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)