

# LH ELISA kit

**Highly sensitive LH ELISA kit enabling detection of LH in serum, plasma and tissue culture media in just 2 hours.**

The LH ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 2 hours.

LH is produced by the gonadotroph cells in the anterior pituitary gland. It is responsible for ovulation in women and controls testosterone synthesis in men. An acute rise of LH, referred to as the “LH surge”, triggers ovulation and development of the corpus luteum in females. After ovulation, LH supports the transient life span of the corpus luteum acting on the luteinized granulosa cells. Cell enlargement and increased progesterone production are involved in the luteinization of the granulosa cells, which are closely associated with the developing gamete (egg) in the ovary.

Citations: 19

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-KIT107-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of LH, detecting as little as 5.2 mIU/ml
- Negligible reactivity from similar glycoprotein hormones
- High throughput format with results in 2 hours for up to 39 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis



## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Luteinizing hormone, Lutropin

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of LH in tissue culture media, serum and plasma in human and rat samples.

**Assay Time** 2 hours

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Microtiter Plate, Assay Buffer 13, Standard, Wash Buffer Concentrate, Antibody, Tracer, Conjugate, Substrate, and Stop Solution

**Crossreactivity** LH (100%), FSH ( $\leq 0.004\%$ ), hCG ( $\leq 0.004\%$ ), TSH ( $\leq 0.3\%$ )

**Sensitivity** 5.2 mIU/ml (1.2-280 mIU/ml)

**Species Reactivity** Human, Rat

**UniProt ID** P01215 ( $\alpha$  chain), P01229 ( $\beta$  chain)

**Wavelength** 450 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)