

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 (α chain), P01229 (β chain)
Wavelength	450 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)