

FSH ELISA kit

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

The FSH ELISA kit is a colorimetric sandwich immunoassay kit with results in 2 hours.

FSH is a glycoprotein produced by the gonadotroph cells in the anterior pituitary gland. It regulates the development, growth, pubertal maturation and reproductive process in both women and men. LH and FSH act synergistically in regulating female ovulation and male spermatogenesis.

Citations: 12

[View Online »](#)

Ordering Information

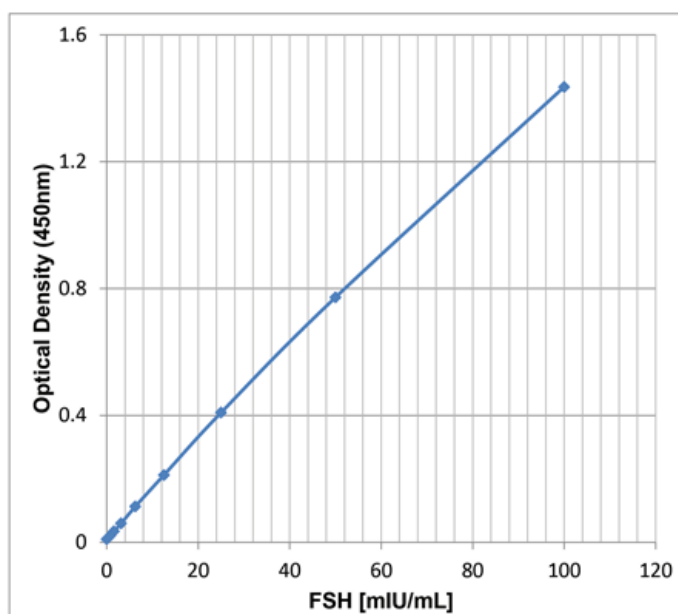
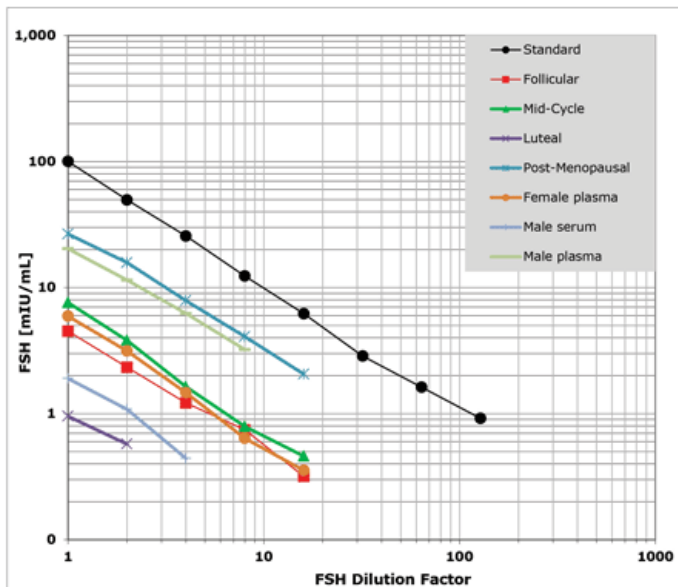
[Order Online »](#)

ENZ-KIT108-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of FSH, detecting as little as 0.5 mIU/ml
- Negligible reactivity from similar glycoprotein hormones
- High throughput format with results in 2 hours for up to 37 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Handling & Storage

Use/Stability	Store all components at +4°C, except the detector antibody, which must be stored at -20°C.
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Follicle-stimulating hormone, Follitropin
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of FSH in tissue culture media, serum and plasma in human 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
Sensitivity	0.5 mIU/ml (0.78-100 mIU/ml)
Species Reactivity	Human
UniProt ID	P01215 (α chain), P01225 (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-
fr@enzolifesciences.com](mailto:info-fr@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)