

ACTH ELISA kit

Highly sensitive ACTH ELISA kit enabling detection of ACTH in human plasma in 4.5 hours.

The ACTH ELISA kit is a colorimetric immunoassay kit with results in 4.5 hours.

Adrenocorticotrophic hormone (ACTH), also known as corticotropin, is a polypeptide tropic hormone produced and secreted by the anterior pituitary gland. It is an important component of the hypothalamic-pituitary-adrenal axis and is often produced in response to biological stress (along with its precursor corticotropin-releasing hormone from the hypothalamus). Its principal effects are increased production and release of corticosteroids

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-KIT138-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of ACTH, detecting as little as 0.46 pg/ml.
- Negligible reactivity to similar hormones.
- High throughput with results in 4.5 hours.
- 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 Adrenocorticotrophic hormone, Corticotropin

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of ACTH in plasma in human samples.

Assay Time 4.5 hours

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

Contents Microtiter Plate, Standard, Wash Buffer Concentrate, Buffer, Antibody, Conjugate, Substrate, and Stop Solution

Sensitivity 0.46 pg/ml (0.5-165 pg/ml)

Species Reactivity Human

UniProt ID P01189 (human)

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)