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.

Adrenocorticotropic 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: 8

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 Adrenocorticotropic hormone, Corticotropin

Application Colorimetric detection, ELISA

Application NotesFor 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

Last modified: May 29, 2024

