

# TSH (human) ELISA kit

Highly sensitive TSH ELISA kit enabling detection of TSH in serum in just 75 minutes.

The TSH ELISA kit is a colorimetric immunometric immunoassay kit with results in 75 minutes.

TSH is a glycoprotein hormone secreted by the anterior pituitary gland and regulates the synthesis and release of  $T_3$  and  $T_4$  by the thyroid gland. TSH has two subunits,  $\alpha$  and  $\beta$ . The  $\beta$  unit gives rise to its biological activity. Serum TSH measurement is one of the most important tools in the diagnosis of thyroid disorders. Increased serum TSH is an early and sensitive indicator of decreased thyroid reserve and overt primary hypothyroidism. Decreased TSH level is an indicator of TSH-independent hyperthyroidism (Graves' disease).

## Ordering Information

[Order Online »](#)

ENZ-KIT131-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of TSH, detecting as little as 0.5  $\mu$ U/ml
- Negligible reactivity from similar glycoprotein hormones
- High throughput format with results in 75 minutes
- 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	Thyroid-stimulating hormone, Thyrotropin, hTSH
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of TSH in serum in human samples.
Assay Time	75 min.
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Microtiter Plate, Standard, Controls, Wash Buffer Concentrate, Conjugate, Substrate, and Stop Solution
Sensitivity	0.5 µIU/ml (0.5-40 µIU/ml)
Species Reactivity	Human
UniProt ID	P01215 (α chain), P01222 (β chain)
Wavelength	450 nm

Last modified: June 25, 2025



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)