

# IL-6 (mouse), ELISA kit

The IL-6 (mouse), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The IL-6 (mouse), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 3 hours.

Interleukin-6 (IL-6) is a cytokine critical to the regulation of the immune and hematopoietic systems. Human IL-6 is expressed by T-cells, mast cells, monocytes, macrophages, fibroblasts, endothelial cells, keratinocytes, and many tumor cell lines. It appears to take part in acute phase reactions and response to injury and inflammation. Human IL-6 also stimulates the differentiation of B-cells for antibody production, promotes expansion of activated T-cells, expands hematopoietic cell production, and induces the expression of acute phase proteins.

Citations: 41

[View Online »](#)

## Ordering Information

[Order Online »](#)

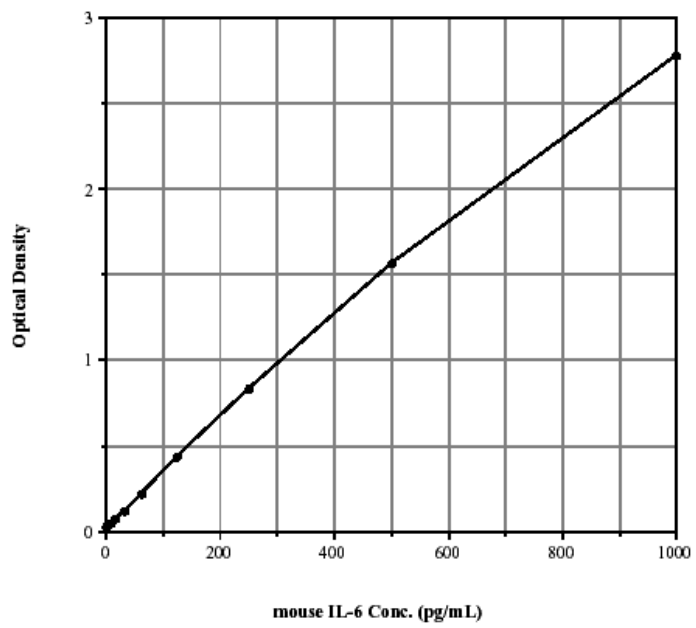
ADI-900-045

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Results in <3 hours
- Sensitive to 1.01 pg/mL
- Measure up to 36 samples per 96 well plate
- Ready-to-use reagents and pre-coated plates reduce errors and save your time



## Handling & Storage

**Use/Stability** Store all components at 4°C, except standard at -20°C.

**Shipping** Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Alternative Name** Interleukin-6

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of mouse Il-6 in culture supernatants and serum. Cited sample type includes tissue.

**Assay Time** <3 hours

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

**Contents** Microtiter plate, Conjugate, Antibody, Assay buffer 13, Wash buffer concentrate, Standard, TMB Substrate, Stop solution

**Sensitivity** 1.01 pg/ml (range 7.81-1,000 pg/ml)

**Species Reactivity** Mouse

**Technical Info / Product Notes** **Cited samples:**  
[Cytokine ELISAs Cited Samples](#)

**UniProt ID** P08505

**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-  
be@enzolifesciences.com](mailto:info-be@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)