

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: 40

[View Online »](#)

Ordering Information

[Order Online »](#)

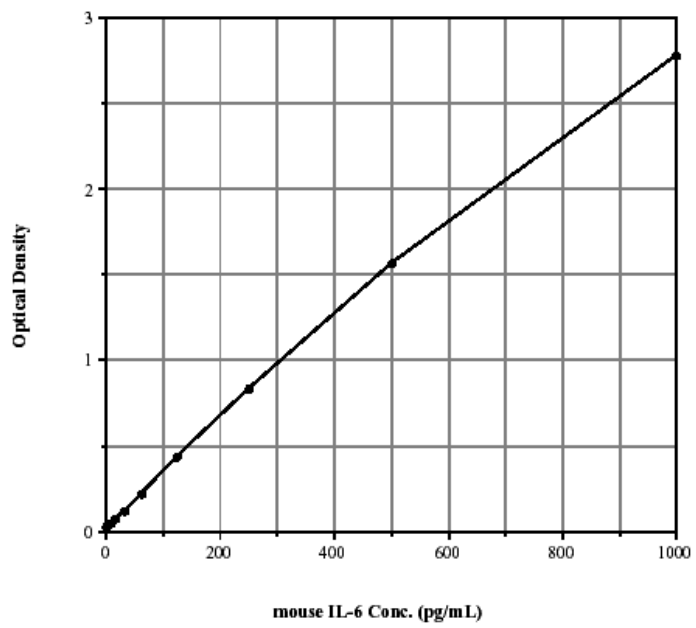
ADI-900-045

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)