

IL-6 (human), high sensitivity ELISA kit

Most sensitive (0.057 pg/mL) colorimetric kit for IL-6 quantification

The IL-6 (human), high sensitivity ELISA kit is a colorimetric immunoassay kit with results in 3 hours. Absorbance is read at 450 nm. Measure up to 40 samples in duplicate per 96 well kit.

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 plays a role 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: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

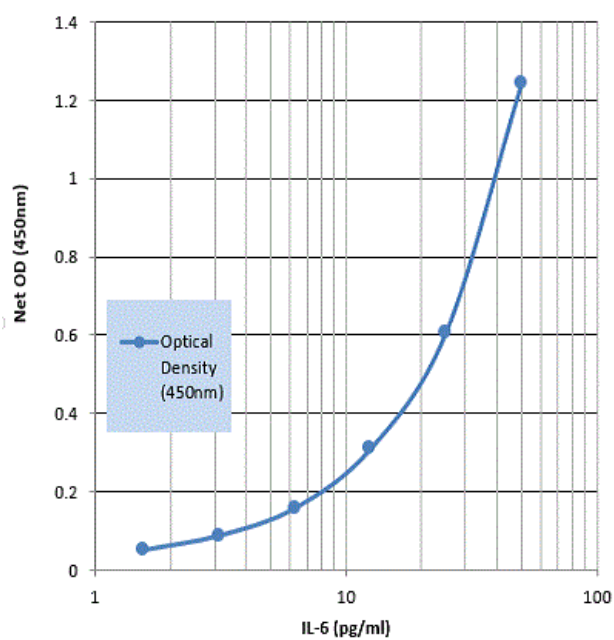
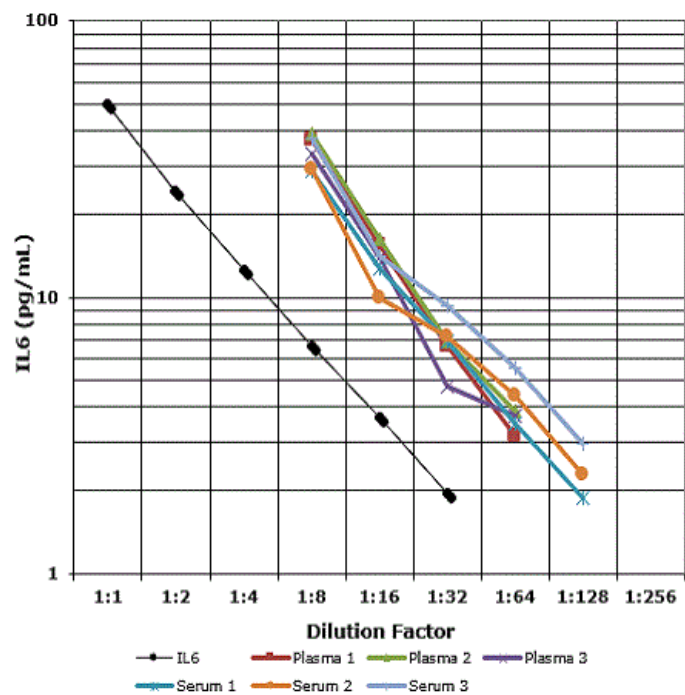
ENZ-KIT178-0001

96 wells

Manuals, SDS & CofA

[View Online »](#)

- High-sensitivity ELISA enables quantification of low amounts of IL-6
- Easy-to-use liquid color-coded reagents reduce errors
- Rapid quantitative results in 3 hours
- High-throughput immunoassay measures up to 40 samples in duplicate per 96 well kit
- Ready-to-use reagents and pre-coated plates reduce errors and save time



Handling & Storage

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

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-6

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human IL-6 in culture supernatants, plasma, serum, and urine.

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

Crossreactivity Negligible crossreactivity to IL-1 α , IL-1 β , IL-2, IL-3, IL-4, IL-7, IL-8, TNF α , and TNF β

Sensitivity 0.057 pg/mL (1.56 – 50 pg/mL)

Species Reactivity Human

UniProt ID P05231

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)