

IL-10 (human), ELISA kit

The IL-10 (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The IL-10 (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 450 nm. Ready-to-use liquid color-coded reagents reduce laboratory errors.

Interleukin-10 (IL-10) was initially identified as the Cytokine Synthesis Inhibitory Factor based on the observation that mouse Th1 cytokine production was suppressed by a Th2-derived factor. IL-10 is a macrophage deactivating factor, acting on macrophage-monocyte accessory cells to produce its inhibitory effects on T cells and natural killer cells. It also regulates growth and/or differentiation of B cells, mast cells, granulocytes, dendritic cells, keratinocytes and endothelial cells. Circulating levels of IL-10 are increased in allergic asthma, systemic sclerosis, a variety of cancers, post-transplantation patients, and in sepsis. The therapeutic potential of IL-10 includes rheumatoid arthritis, systemic lupus erythematosus, multiple sclerosis, and HIV infection.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

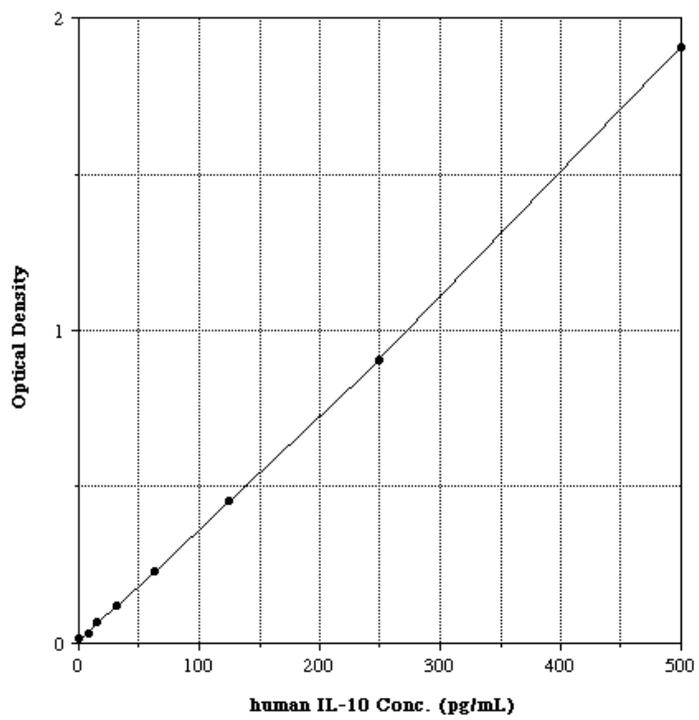
ADI-900-036

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use - liquid color-coded reagents reduce error
- Sensitive - detect < 4 pg/ml of IL-10
- Fast - results in < 3 hours from up to 39 samples in duplicate
- 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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-10

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human IL-10 in culture supernatants, plasma, and serum.

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 3.75 pg/ml (range 7.81 – 500 pg/ml)

Species Reactivity Human

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

UniProt ID P22301

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-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)