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: 8

View Online »

Ordering Information

Order Online »

ADI-901-036

5x96 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 precoated plates reduce errors and save your time
- Save with bulk kit package!

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.

Assay Time <3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents Microtiter plate, Conjugate, Antibody, Assay buffer 13

concentrate, Wash buffer concentrate, Standard, TMB

Substrate, Stop solution 2

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com