

IL-1 β (human), ELISA kit

The IL-1 β (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The IL-1 β (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 450 nm. Quantify 1 pg/mL IL-1 β with lot sample volume needed saving you precious sample.

The interleukin-1 (IL-1) cytokine family consists of IL-1 α , IL-1 β , and IL-1R α (receptor antagonist) which are expressed primarily by endothelial cells and activated macrophages. IL-1 α and IL-1 β elicit pro-inflammatory cytokine production and cellular activation, and are the primary mediators of hypothalamic regulation that results in fever. IL-1 α and IL-1 β signals are transmitted via a dual subunit IL-1 receptor composed of IL-1R1 and the IL-1 receptor accessory protein (IL-1RAcP). IL-1RAcP is vital to IL-1 signaling, as it recruits the IL-1 receptor-associated kinase (IRAK) to the receptor complex, initiating NF- κ B pathway activation.

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-900-130A

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive - quantify down to 1 pg/mL
- Efficient - only 50 μ l diluted sample required saving precious materials
- Specific - no cross-reactivity with related cytokines
- Results from up to 40 samples in duplicate in just 3 hours
- Ready-to-use reagents and pre-coated plates reduce errors and save your time

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-1 β

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of IL-1 β in culture supernatant, plasma, serum, and urine of human origin.

Assay Time < 3 hours

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

Contents Standard, Standard Diluent Buffer, Coated 96-well plate, Conjugate, Streptavidin-HRP, Streptavidin-HRP Diluent, Wash Buffer Concentrate, TMB Solution, Stop Solution, Plate Sealers

Crossreactivity The following substances were tested and found to have no cross-reactivity: human IL-1 α , IL-1ra, IL-2, IL-3, IL-4, IL-6, IL-7, IL-8, IL-10, IFN- α , TGF- β , GM-CSF, G-CSF, OSM, MIP-1 α , MIP-1 β , LIF, MCP-1, G-CSF, RANTES, PF-4, β TG, GRO, IP-10, SCF, TNF- α , and TNF- β .

Sensitivity 1 pg/mL

Species Reactivity Human

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

UniProt ID P01584



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

450 nm

European Sales Office
ENZO LIFE SCIENCES
ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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