

# IL-1 $\beta$ (rat), ELISA kit

The IL-1 $\beta$  (rat), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The IL-1 $\beta$  (rat), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 4 hours.

The interleukin-1 (IL-1) cytokine family consists of IL-1 $\alpha$ , IL-1 $\beta$ , and IL-1Ra (receptor antagonist) which are expressed primarily by endothelial cells and activated macrophages. IL-1 $\alpha$  and IL-1 $\beta$  elicit pro-inflammatory cytokine production and cellular activation, and are the primary mediators of hypothalamic regulation that results in fever. IL-1 $\alpha$  and IL-1 $\beta$  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- $\kappa$ B pathway activation.

Citations: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-900-131A

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Sensitive - use less precious sample per test
- Higher throughput - test up to 40 samples in duplicate in less than 4 hours
- Quantitative - obtain statistically significant results and compare data between multiple experiments
- Ready-to-use - pre-coated plates save time and reduce errors

# Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

# Regulatory Status

 RUO - Research Use Only

# Product Details

Alternative Name

Interleukin-1 $\beta$

Application

Colorimetric detection, ELISA

Application Notes

For the quantitative determination of IL-1 $\beta$  in culture supernatants and serum of rat origin.

Assay Time

<4 hours

Compatibility

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

Contents

Microtiter plate, SA-HRP concentrate, SA-HRP dilution buffer, Biotinylated antibody, Standard, Standard diluent, Wash buffer concentrate, TMB Substrate, Stop solution

Sensitivity

4 pg/mL (range 31.3 – 2,000 pg/mL)

Species Reactivity

Rat

Technical Info / Product Notes

**Cited samples:**  
[Cytokine ELISAs Cited Samples](#)

UniProt ID

Q63264

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)