

IL-1 β (rat), ELISA kit

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

The IL-1 β (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 α , IL-1 β , and IL-1Ra (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: 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 β
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of IL-1 β 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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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