

# IL-2 (human), ELISA kit

The IL-2 (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The IL-2 (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours.

Interleukin-2 (IL-2) plays a critical role in the mediation of immune response by acting as a growth and differentiation factor for T cells, as well as B cells, natural killer (NK) cells, lymphokine-activated killer (LAK) cells, tumor-infiltrating lymphocytes (TIL), monocytes, macrophages, and oligodendrocytes. The human IL-2 receptor can be composed of several combinations of a 55 kDa IL-2R $\alpha$  subunit, a 70-75 kDa IL-2R $\beta$  subunit, and a 64 kDa IL-2R $\gamma$  subunit. IL-2 binding affinity varies depending on which combination of subunits is present in the bonding complex, with the high affinity binding site having all three subunits. Studies have shown IL-2 to cause the regression of some types of tumors and it may have implications in immunotherapy and the treatment of infectious diseases and immune deficiencies.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-900-118A

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Results in just 3 hours from up to 39 samples in duplicate
- Low dynamic range provides sensitive measurements
- Validated in a wide variety of matrices
- 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** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Interleukin-2

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of human IL-2 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, Wash buffer concentrate, Standard, TMB Substrate, Stop solution 2

**Sensitivity** 6.6 pg/ml (range 7.81 – 500 pg/ml)

**Species Reactivity** Human

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

**UniProt ID** P60568

**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-  
be@enzolifesciences.com](mailto:info-be@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-  
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)