

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 α subunit, a 70-75 kDa IL-2R β subunit, and a 64 kDa IL-2R γ 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.

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)