

IL-2 (human), (recombinant)

Interleukin 2 (IL-2) is an immunomodulatory cytokine that is produced by lymphocytes. IL-2 signals through the IL-2R receptor to induce activated T cell proliferation and promote T cell differentiation. IL-2 also stimulates the proliferation and differentiation of B cells, natural killer cells, monocytes, and macrophages.

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-PRT254-0100 | 100µg |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)

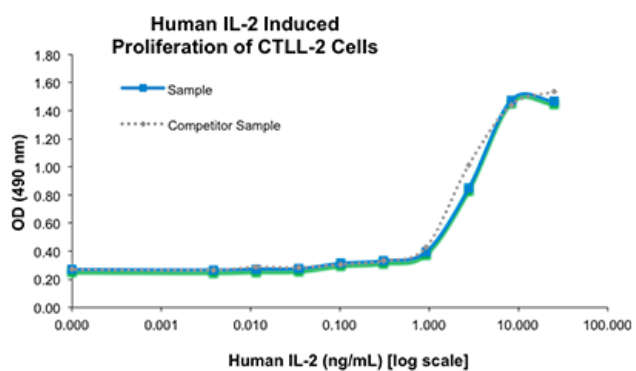


Figure 1: CTLL-2 cells were cultured with 0 to 25 ng/mL Human IL-2. Cell proliferation was measured after 66 hours and the linear portion of the curve can be used to calculate the ED50.

Handling & Storage

| | |
|-------------------|---|
| Handling | Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution. |
| Long Term Storage | -80°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------|---|
| Alternative Name | Interleukin-2, T cell growth factor, TCGF, Aldesleukin |
| Endotoxin Content | ≤1 EU/μg protein (LAL test) |
| Formulation | Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% Trifluoroacetic Acid (TFA). |
| MW | ~15.5kDa |
| Purity | ≥95% (SDS-PAGE) |
| Reconstitution | Reconstitute in 10 mM HCl at 0.1 mg/mL. |
| Source | Produced in <i>E. coli</i> . Contains 134 amino acids. |
| UniProt ID | P60568 |

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)