

IL-15 (human), (recombinant)

Interleukin 15 (IL-15) is a widely expressed proinflammatory cytokine that is structurally and functionally related to interleukin 2 (IL-2). IL-15 signals through JAK kinases to activate STAT3, STAT5, and STAT6 transcription factors. IL-15 regulates the activation of T cells, neutrophils, and macrophages, and is a stimulatory cytokine promoting dendritic cell function. IL-15 expression is dysregulated in chronic inflammatory disorders, including rheumatoid arthritis, psoriasis, and pulmonary inflammatory diseases.

Ordering Information

[Order Online »](#)

ENZ-PRT255-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Human IL-15 Induced Proliferation of CTLL-2 Cells

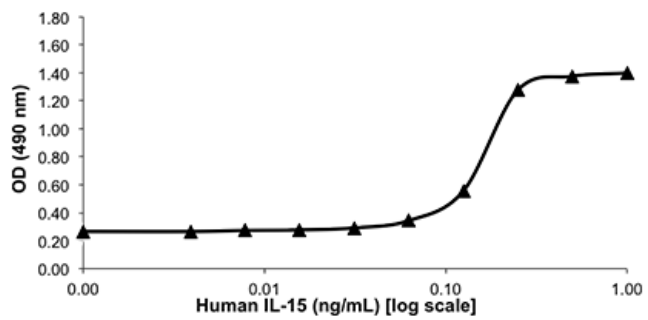


Figure 1: Serial dilutions of Human IL-15 (starting at 1 ng/mL) were added to CTLL-2 cells with other needed cytokines and stimuli. After 65 hours, cell viability was measured and the linear portion of the curve can be used to calculate the ED50.

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-15, IL-T
Endotoxin Content	≤1 EU/μg protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 10 mM sodium bicarbonate, pH 8.
MW	~12.9kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 115 amino acids.
UniProt ID	P40933



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)