

IL-21 (human), (recombinant)

Interleukin 21 (IL-21) is normally expressed by activated CD4+ T cells. The IL-21 receptor activates the JAK/STAT signaling pathway and is expressed on T, B, and natural killer (NK) cells. IL-21 cooperates with IL-4 for the production of multiple antibody classes in B cells. IL-21 has pleiotropic effects on the proliferation, differentiation, and effector functions of B, T, NK, and dendritic cells.

Ordering Information

[Order Online »](#)

ENZ-PRT248-0010	10µg
ENZ-PRT248-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

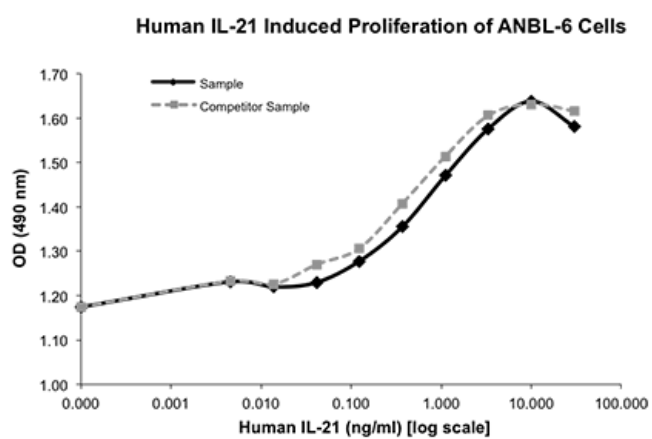


Figure 1: Serial dilutions of Human IL-21 (starting at 30 ng/mL) were added to ANBL-6 cells. After 84 hours, cell proliferation was measured and the linear portion of the curve can be used to calculate the ED50.

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-21
Biological Activity	B9 cell proliferation; ED50 ≤ 50 ng/mL ($\geq 2.0 \times 10^4$ units/mg).
Endotoxin Content	≤1 EU/μg protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 10 mM sodium phosphate, pH 7.5.
MW	~15.4 kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 132 amino acids.
UniProt ID	Q9HBE4



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)