

IL-22 (human), (recombinant)

Interleukin 22 (IL-22), also called IL-TIF, is an IL-10 family member that is produced by activated dendritic cells and T lymphocytes. IL-22 signals via a heteroduplex receptor consisting of IL-22R and IL-10RB chains. IL-22 is a potent mediator of cellular inflammatory responses.

Ordering Information

[Order Online »](#)

ENZ-PRT250-0010	10µg
ENZ-PRT250-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

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-22, IL-TIF, IL-10-related T-cell-derived-inducible factor, Cytokine Zcyto18
Endotoxin Content	≤1 EU/μg protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 10 mM sodium citrate, pH 3.0.
MW	Dimer, ~16.9/33.8 kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 147/294 amino acids.
UniProt ID	Q9GZX6



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com