

# IL-22 (mouse), (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.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

|                          |   |
|--------------------------|---|
| <b>Handling</b>          | 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. |
| <b>Long Term Storage</b> | -80°C   |
| <b>Shipping</b>          | Ambient Temperature   |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                            |  |
|----------------------------|--|
| <b>Alternative Name</b>    | Interleukin-22, IL-TIF, IL-10-related T-cell-derived-inducible factor, Cytokine Zcyto18                |
| <b>Biological Activity</b> | Production of IL-10 from human COLO-205 cells; ED50 $\leq 1$ ng/mL ( $\geq 1.0 \times 10^6$ units/mg). |
| <b>Endotoxin Content</b>   | $\leq 1$ EU/ $\mu$ g protein (LAL test)  |
| <b>Formulation</b>         | Lyophilized from a sterile (0.2 micron) filtered solution containing 0.1% Trifluoroacetic Acid (TFA).  |
| <b>MW</b>                  | ~16.8 kDa  |
| <b>Purity</b>              | $\geq 95\%$ (SDS-PAGE)   |
| <b>Reconstitution</b>      | Reconstitute in sterile water at 0.1 mg/mL.  |
| <b>Source</b>              | Produced in <i>E. coli</i> . Contains 147 amino acids.   |
| <b>UniProt ID</b>          | Q9JJY9   |

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)