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 *E. coli*. Contains 147/294 amino acids.

UniProt ID Q9GZX6

Last modified: May 29, 2024

