

I882-40N-10	10 µg
I882-40N-100	100 µg
I882-40N-1000	1 mg

Recombinant Human IL22

Recombinant human protein expressed in E. coli cells

Catalog # I882-40N

Lot # G3211-68

Product Description

Recombinant Human IL22 was expressed in E. coli cells. The protein accession number is [Q9GZX6](#)

Gene Aliases

IL-22, IL-TIF, Cytokine Zcyto18, IL-10-related T cell-derived-inducible factor, ILTIF, ZCYTO18

Recombinant Human Interleukin-22 (IL22)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

Formulation

Recombinant Interleukin-22 was lyophilized from 0.2 µm filtered concentrated (1.0 mg/mL) solution in 500 mM NaCl, 20 mM Tris, pH 8.0.

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

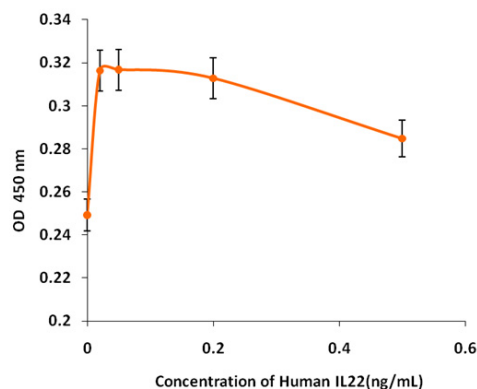
Storage and Stability

Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

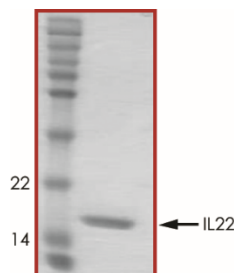
Originally identified as a gene induced by IL-9 in mouse T cells and mast cells, IL-22 is a member of the IL-10 family of cytokines. IL-22 inhibits IL-4 production by Th2 cells, and induces acute phase reactants in the pancreas and liver. The IL-22 receptor complex consists of two receptor subunits, IL-22R and IL-10R-beta that belong to the class II cytokine receptor family. Recombinant human IL-22 is a non-disulfide-linked homodimeric protein.

Specific Activity



Activity was determined by the ability to inhibit growth of human A549 cancer cells at concentrations ranging between 50-100 ng/mL.

Purity



Recombinant Human IL22 resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.0 kDa**

Recombinant Human IL22

Recombinant human protein expressed in E. coli cells

Catalog Number	I882-40N
Specific Lot Number	G3211-68
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.