

Recombinant Human IL13

Recombinant human protein expressed in E. coli cells

Catalog # **1873-40N**

Lot # **G3211-56**

Product Description

Recombinant Human IL13 was expressed in E. coli cells. The protein accession number is [P35225](#)

Gene Aliases

IL-13, NC30

Recombinant Human Interleukin-13 (IL13)

Endotoxin Level

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

Formulation

Lyophilized from a 0.2 µm filtered solution in sodium phosphate and NaCl (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.

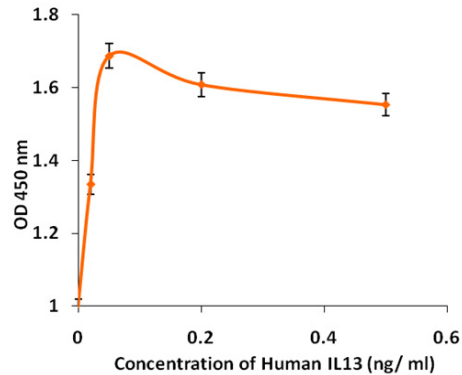
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. 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

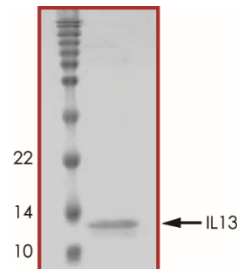
Human IL13 is expressed in activated T helper cells (resembling Th0, Th1, Th2), and T cells expressing CD8. The alpha chain of the IL13 receptor has weak binding activity for IL13. High affinity receptors are formed when the receptor alpha chains of IL13 and IL4 receptors are coexpressed. Different IL13 receptor structures have been shown to exist on various cell types and the IL13 receptor may share more than one component with IL4 receptor. IL13 down-modulates macrophage activity, reducing the production of pro-inflammatory cytokines and chemokines in response to IFN-gamma or bacterial lipopolysaccharides. IL13 enhances the production of the IL1 receptor antagonist IL1ra. IL13 also decreases the production of nitric oxide by activated macrophages, leading to a decrease in parasitocidal activity. IL13 induces differentiation of human monocytes, enhances survival time in culture, and also induces differentiation and proliferation and isotype switching in B cells.

Specific Activity



The ED50, as determined by the dose-dependent proliferation of TF-1 cells was found to be ≤ 2.0 ng/mL

Purity



Recombinant Human IL13 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **13.0 kDa**

Recombinant Human IL13

Recombinant human protein expressed in E. coli cells

Catalog Number	1873-40N
Specific Lot Number	G3211-56
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. 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.