

Recombinant Human IL3

Recombinant human protein expressed in *E. coli* cells

Catalog # 1863-40N

Lot # G3211-70

Product Description

Recombinant Human IL3 was expressed in *E. coli* cells. The protein accession number is [P08700](#)

Gene Aliases

IL-3, Mast cell growth factor, MCGF, Multi-CSF, Hematopoietic growth factor, HCGF, P-cell stimulation factor
Recombinant Human Interleukin-3 (IL3)

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 Phosphate Buffer and NaCl (pH 6.3)

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

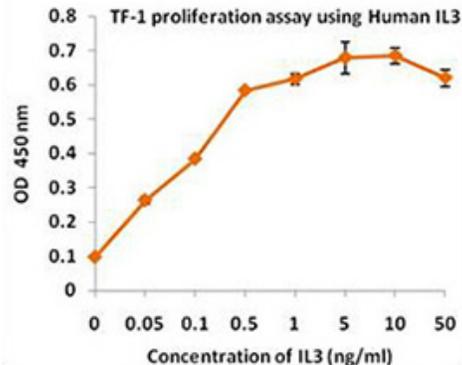
Storage and Stability

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

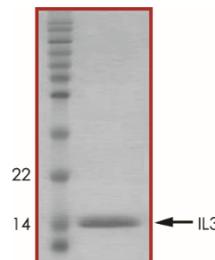
IL3 is produced mainly by T cells following cell activation by antigens and mitogens, but also by keratinocytes, natural killer cells, mast cells, endothelial cells, and monocytes. The analysis of bacterial- derived recombinant IL3 shows that glycosylation is not required for activity. IL3 sequences are evolutionarily less well conserved with human and murine IL3 sharing approximately 29% homology (at the protein level) and murine and rat IL3 sharing approximately 54% homology. IL3 receptors are expressed on macrophages, mast cells, eosinophils, megakaryocytes, basophils, bone marrow progenitor cells and various myeloid leukemia cells. Binding of IL3 to its receptor causes specific phosphorylation of a 150 kDa membrane glycoprotein. Recombinant human IL3 is a non-glycosylated globular protein.

Specific Activity



The ED50 as determined by the dose-dependent proliferation of Human TF-1 cells was found less than 0.5 ng/mL

Purity



Recombinant Human IL3 resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **15.0 kDa**

Recombinant Human IL3

Recombinant human protein expressed in *E. coli* cells

Catalog Number	1863-40N
Specific Lot Number	G3211-70
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.