

Recombinant Human IL5

Recombinant human protein expressed in E. coli cells

Catalog # 1865-40N

Lot # G3211-77

Product Description

Recombinant Human IL5 was expressed in E. coli cells. The protein accession number is [P05113](#)

Gene Aliases

IL-5, B cell growth factor II, BCGF-II, Cytotoxic T lymphocyte inducer, Eosinophil differentiation factor T cell replacing factor, TRF Recombinant Human Interleukin-5 (IL5)

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 Tris with NaCl

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.

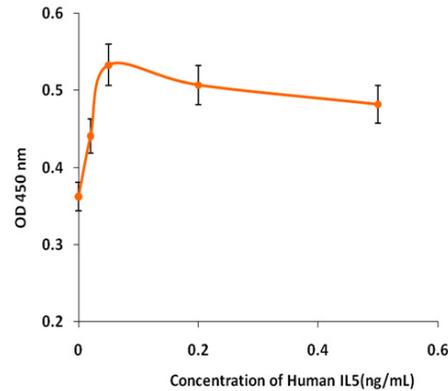
Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favourable performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

Interleukin 5 (IL5) is a specific hematopoietic growth factor that is responsible for the growth and differentiation of eosinophils. IL5 promotes the growth of immature hematopoietic progenitor cells BFU-E while it causes differentiation of CFU-E the proliferation of which is inhibited by IL5. B cells can be made responsive to IL5 by treatment with suboptimal doses of IL1. Interleukin 5 also promotes the generation of cytotoxic T cells from thymocytes. In thymocytes IL5 induces the expression of high affinity IL2 receptors. Human and mouse IL5 is cross-species reactive. Recombinant Human IL5 is a 26.5 kDa disulfide-linked homodimerized protein.

Specific Activity



Purity



Recombinant Human IL5 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-251

Approx. MW **13.0 kDa**

Recombinant Human IL5

Recombinant human protein expressed in E. coli cells

Catalog Number	1865-40N
Specific Lot Number	G3211-77
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favourable performance, avoid repeated handling and multiple 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.