

Recombinant Human IL1B

Recombinant human protein expressed in E. coli cells

Catalog # I861-40BN

Lot # G3211-65

Product Description

Recombinant Human IL1B was expressed in E. coli cells. The protein accession number is [P01584](#)

Gene Aliases

Catabolin, Lymphocyte-activating factor, LAF, Endogenous Pyrogen, EP, IL1F2, Recombinant Human Interleukin-1 Beta (IL1B)

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, NaCl, EDTA and DTT (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.

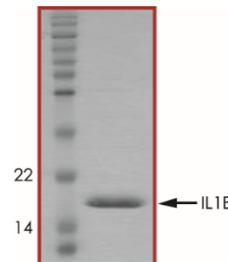
Scientific Background

Monocytes are the main source of secreted IL1; expressing mostly IL1B, while human keratinocytes express large amounts of IL1A. Overall, IL1B is the predominant form in humans while, IL1A is the one in mice. Both isoforms of IL1 bind to the same receptor and therefore show similar, if not identical, biological activities. The IL1B precursor though must be processed before it can bind to the receptor. IL1 proteins are involved in the inflammatory response, being identified as endogenous pyrogens, and are reported to stimulate the release of prostaglandin and collagenase from synovial cells. A monomer, IL1B stimulates thymocyte proliferation, B-cell maturation and proliferation and fibroblast growth factor activity.

Specific Activity

The ED50 as determined by the dose-dependent inhibition of prostate carcinoma cell line is less than 1ng/ml

Purity



Recombinant Human IL1B resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **17.5 kDa**

Recombinant Human IL1B

Recombinant human protein expressed in E. coli cells

Catalog Number	I861-40BN
Specific Lot Number	G3211-65
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.