

## Recombinant Human IL1A

Recombinant human protein expressed in E. coli cells

**Catalog # I861-40N**

Lot # G3211-64

### Product Description

Recombinant Human IL1A was expressed in E. coli cells. The protein accession number is [P01583](#)

### Gene Aliases

Hematopoietin-1, Lymphocyte-activating factor, LAF, Mononuclear Cell Factor, MCF, IL1F1, Recombinant Human Interleukin-1 Alpha (IL1A)

### 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 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.

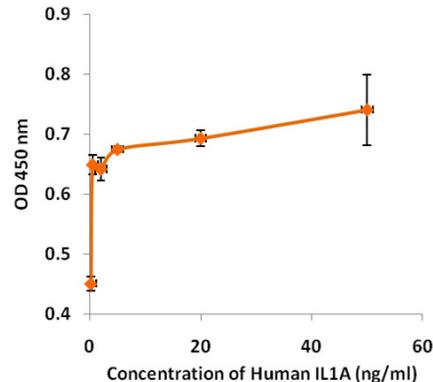
### Storage and Stability

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

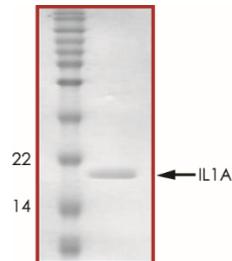
There are two functionally comparable forms of IL1: IL1A and IL1B, that are encoded by two different genes. IL1B is the predominant form in humans while, it is IL1A in mice. Both forms of IL1 bind to the same receptor and therefore show similar, if not identical, biological activities. Latter include but not limited to, stimulation of T-helper cells, promotion, proliferation and synthesis of immunoglobulins and the activation of fibroblasts and NK cells. Unlike IL1B, the IL1A precursor does not require processing before it can bind to the receptor. Monocytes are the main source of secreted IL1B while human keratinocytes are known to express large amounts of IL1A. Recombinant Human IL1A is a non-glycosylated polypeptide of 18 kDa.

### Specific Activity



The ED50 as determined by the dose-dependent stimulation of murine D10S cells was found to be ≤ 0.002 ng/mL

### Purity



Recombinant Human IL1A resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **18.0 kDa**

## Recombinant Human IL1A

Recombinant human protein expressed in E. coli cells

Catalog Number	I861-40N
Specific Lot Number	G3211-64
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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