

| Catalog #    | Aliquot Size |
|--------------|--------------|
| I866-40N-20  | 20 µg        |
| I866-40N-100 | 100 µg       |

## Recombinant Human IL6

Recombinant human protein expressed in E. coli cells

**Catalog # I866-40N**

Lot # G3211-78

### Product Description

Recombinant Human IL6 was expressed in E. coli cells. The protein accession number is [P05231](#)

### Gene Aliases

IL-6, B cell differentiation factor, BCDF, B cell stimulatory factor 2, BSF-2, CTL differentiation factor, CDF, HPGF, HSF, MGI-2, Hybridoma growth factor, Interferon beta-2, IFN-beta 2, Recombinant Human Interleukin-6 (IL6)

### 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 PBS (pH 7.4)

### Reconstitution Protocol

A quick spin of the vial followed by reconstitution in sterile PBS to a concentration not less than 0.1 mg/mL.

### 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.

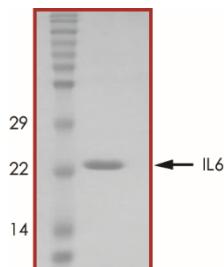
### Scientific Background

IL6 is a pleiotropic cytokine that participates in a wide variety of biological functions such as acute phase response, inflammation, hematopoiesis, bone metabolism and cancer. While produced mainly by stimulated monocytes, fibroblasts and endothelial cells, IL6 is also known to be produced by macrophages, T cells, B lymphocytes, granulocytes, smooth muscle cells, eosinophils, chondrocytes, osteoblasts, mast cells, glial cells and keratinocytes, upon stimulation. Unlike murine IL6 which is inactive on human cells, both human and murine IL6 are equally active on murine cells. The IL6 receptor is a strongly glycosylated protein of 80 kDa and a length of 449 amino acids (designated CD126). Recombinant human IL6 is a 21 kDa, non-glycosylated protein that contains two disulfide bridges.

### Specific Activity

The ED50 as determined by the dose-dependent stimulation of human TF-1 cells was in the range of 0.5-1 ng/mL

### Purity



Recombinant Human IL6 resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **21.0 kDa**

## Recombinant Human IL6

Recombinant human protein expressed in E. coli cells

|                     |   |
|---------------------|---|
| Catalog Number      | I866-40N  |
| Specific Lot Number | G3211-78  |
| 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.**