

## Recombinant Human IP-10 (CXCL10)

Recombinant human protein expressed in E. coli cells

**Catalog # 1838-40BN**

Lot # G3211-83

### Product Description

Recombinant Human IP-10 (CXCL10) was expressed in E. coli cells. The protein accession number is [P02778](#)

### Gene Aliases

CXCL10, C-X-C motif chemokine 10, 10 kDa interferon gamma-induced protein, Gamma-IP10, Small-inducible cytokine B10, IP10

### Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method.

### Formulation

Lyophilized from 0.2 µm filtered PBS solution, pH 7.0.

### 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.

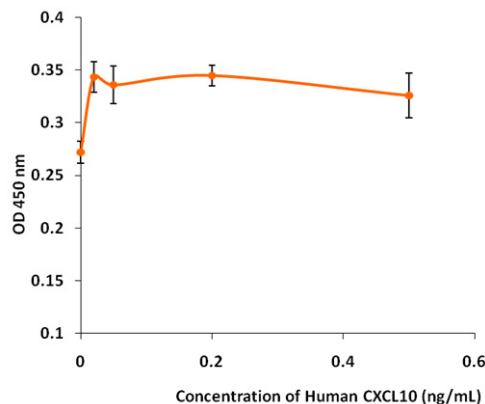
### Storage and Stability

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

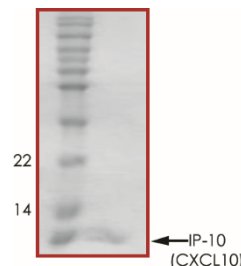
IP-10 shows homology to PF-4 (platelet factor-4) and belongs to the family of chemotactic cytokines known as chemokines. The receptor for IP-10 is CXCR3. IP-10 has been shown to bind to the virus-encoded viroreceptor M3. The expression of IP-10 from a variety of cells, including monocytes, endothelial cells, keratinocytes, and fibroblasts, is induced by IFN-gamma and TNF-alpha. Human neutrophils produce IP-10 in response to IFN-gamma in combination with either TNF-alpha or bacterial lipopolysaccharides. IP-10 probably also plays a role in regulation of the growth of immature hematopoietic progenitor cells. It has been shown to suppress in vitro colony formation of highly enriched cells expressing the cell surface marker CD34 in the presence of SCF, GM-CSF, or SCF and EPO, but not in their absence with the exception of SCF.

### Specific Activity



Determined by its ability to chemoattract human T-Lymphocytes using a concentration of 10-50 ng/mL.

### Purity



Recombinant Human IP10 (CXCL10) resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **9.0 kDa**

## Recombinant Human IP-10 (CXCL10)

Recombinant human protein expressed in E. coli cells

Catalog Number	1838-40BN
Specific Lot Number	G3211-83
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.**