

**Description:** Recombinant Interferon gamma is a disulfide-linked monomer protein consisting of 135 amino acid residue subunits, and migrates as an approximately 16 kDa protein under non-reducing and reducing conditions.

**Source:** Optimized DNA sequence encoding rat Interferon gamma mature chain expressed in E. coli.

**Formulation:** Lyophilized from a 0.2 µm filtered PBS solution, pH 7.0.

**Reconstitution:** 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:** The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8° C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

**Purity:** >95%, as determined by SDS-PAGE and HPLC.

**Endotoxin Level:** Endotoxin level was found to be < 0.1 ng/µg (1EU/µg), using the LAL gel clot method.

**Biological Activity:** The ED(50) was determined by the cytopathic inhibition assay with murine L929 cells infected by EMC virus, and was found to be < 1 ng/ml, corresponding to a specific activity of > 1 x 10<sup>6</sup> units/mg.

**Protein Sequence:**

MSATRRVLVLQLCLMALSGC  
YCQGLTIESLESLSKKNYFNSS  
SMDAMEGKSLLLDIWRNWQK  
DGNTKILESQIISFYLRIFE  
VLKDNQAISNNISVIESHLI  
TNFFSNSKAKKDAFMSIAKF  
EVNNPQIQHKAVNELIRVIH  
QLSPESSLRKRKRSRC

**References:**

1. Sikorski K, et al. *Cytokine Growth Factor Rev.* 2011 Aug,22(4):211-9.
2. Gysemans C, et al. *Biochem Soc Trans.* 2008 Jun,36(Pt 3):328-33.
3. Rameshwar P. *Cell Res.* 2008 Aug,18(8):805-6.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)