

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 10⁶ units/mg.

Protein Sequence:

MSATTRVLVLQLCLMALSGC
YCQGLIESLESLSKKNYFNSS
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

Please visit our website at: www.bpsbioscience.com