



6405 Mira Mesa Blvd, Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Human Interferon-alpha 2b Recombinant

Catalog #: 90159-B

Lot#: 220718

Description: Recombinant Interferon- α 2b is a disulfide-linked monomer protein consisting of 166 amino acid residues, migrating as an approximately 19 kDa protein under non-reducing and reducing conditions.

Source: Optimized DNA sequence encoding Interferon- α 2b mature chain was expressed in *E. coli* cells.

Formulation: Lyophilized from a 0.2 μ m-filtered solution of PBS, 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: >98%, as determined by SDS-PAGE and HPLC.

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

Biological Activity: Determined by a viral resistance assay by to be $\geq 5 \times 10^8$ IU/mg.

Protein Sequence:

MALTFYLLVALVLSYKSFSSLGCDLQTHSLGSRRTLMLLAQMRKISLFSCLKDRHDFG
FPQEEFGNQFQKAETIPVLHEMIQQIFNLFSTKDSSAAWDETLLDKFYTELYQQLNLDL
ACVIQGVGVTEPLMKEDSILAVRKYFQRITLYLKEKKYSPCAWEVVRAEIMRSFSLSTN
LQESLRSKE

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