



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 2a Recombinant

Catalog #: 90158-A
Lot#: 230818

Description: Recombinant Interferon- α 2a 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 2a 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 $2 - 8^{\circ}\text{C}$ 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 $< 0.1 \text{ ng}/\mu\text{g}$ (1EU/ μg), using the LAL gel clot method.

Biological Activity: Determined by a viral resistance assay of human WISH cells to be in the range of 1×10^8 IU/mg.

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