



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

## Human Interferon-alpha 2a Recombinant

Catalog #: 90158-B  
Lot#: 230818

**Description:** Recombinant Interferon- $\alpha$  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- $\alpha$  2a mature chain was expressed in *E. coli* cells.

**Formulation:** Lyophilized from a 0.2  $\mu$ 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  $-20^{\circ}\text{C}$ . Upon reconstitution, store in working aliquots at  $2 - 8^{\circ}\text{C}$  for up to one month, or at  $-20^{\circ}\text{C}$  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/ $\mu\text{g}$ ), using the LAL gel clot method.

**Biological Activity:** Determined by a viral resistance assay of human WISH cells to be in the range of  $1 \times 10^8$  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](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)