

# IFN $\beta$ 1a

Recombinant human interferon beta 1 a

Catalog # rhifn-b1a

For research use only-not for diagnostic or therapeutic use

Version # 10B19-MM

## PRODUCT INFORMATION

### Content:

- IFN $\beta$ 1a is provided as 10  $\mu$ g of a sterile filtered white lyophilized powder.

### Storage and stability:

- IFN $\beta$ 1a is shipped at room temperature. Upon receipt it should be stored at -20°C although it is stable at room temperature for 3 weeks.

- Upon resuspension, IFN $\beta$ 1a should be aliquoted and stored at 4°C for short term storage or -20°C for long term storage. Product is stable one week at 4°C and 12 months at -20°C when properly stored. For long term storage, it is recommended to add a carrier protein (0.1% HSA or BSA).

*Note: Avoid freeze-thaw cycles.*

## DESCRIPTION

Recombinant human IFN-beta 1a (IFN $\beta$ 1a) produced in Chinese Hamster Ovarian (CHO) cells is a single, glycosylated, polypeptide chain containing 166 amino acids and having a molecular mass of 22500 Daltons. Recombinant human IFN $\beta$ 1a is purified by proprietary chromatographic techniques.

### Quality control

- Purity greater than 98% as determined by:

- (a) Analysis by RP-HPLC
- (b) Analysis by SDS-PAGE

- Protein quantification was carried out by two independent methods:

1. UV spectroscopy at 280 nm.
2. Analysis by HLPC, using a calibrated solution of IFN-beta as a reference Standard.

## CHARACTERISTICS

**Amino Acid Sequence:** The sequence of the first five N-terminal amino acids was determined and was found to be Ser-Tyr-Asn-Leu-Leu.

**Specific Activity:** 270 x 10<sup>6</sup> IU/mg as determined in a viral resistance assay using human "Wish" cell line and VSV virus or the monkey VERO cell line with EMCV virus.

**Working Concentration:** 10- 10<sup>4</sup> IU/ml.

**Endotoxin:** Less than 0.1 ng/ $\mu$ g (IEU/ $\mu$ g) of IFN $\beta$ 1a.

**Formulation:** IFN $\beta$ 1a was lyophilized from a 1 mg/ml PBS solution (pH 7.4).

**Solubility:** It is recommended to reconstitute lyophilized IFN $\beta$ 1a in sterile water not less than 100  $\mu$ g/ml, which can then be further diluted to other aqueous solutions.

---

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873  
Outside US: (+1) 858-457-5873  
Europe: +33 562-71-69-39  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)  
Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite A  
San Diego, CA 92121 - USA