

# Recombinant human IFN- $\alpha$ 10

Mammalian cell-expressed human interferon alpha 10 (alpha C) with HSA

Cat. code: rcyc-hifna10

<https://www.invivogen.com/human-ifna2b>

For research use only

Version 23A16-NJ

## PRODUCT INFORMATION

### Contents

- 1  $\mu$ g of lyophilized recombinant human IFN- $\alpha$ 10

Note: This product is sterile filtered prior to lyophilization.

- 1.5 ml endotoxin-free water

### Storage and stability

• Recombinant human IFN- $\alpha$ 10 is shipped at room temperature. Upon receipt it should be immediately stored at -20 °C. Lyophilized product is stable for 6 months at -20 °C.

- Upon resuspension, prepare aliquots of recombinant human IFN- $\alpha$ 1 and store at -80 °C for 4 months.

Note: Avoid repeated freeze-thaw cycles.

### Quality control

- Purity:  $\geq$  95% (SDS-PAGE)
- Endotoxin:  $\leq$  1 EU/ $\mu$ g
- The biological activity has been confirmed using HEK-Blue™ IFN- $\alpha$ / $\beta$  cells (see validation data sheet available on our website).
- The units have been determined for each lot using a corresponding calibrated standard.

### Characteristics

- Source: Mammalian; Chinese hamster ovary (CHO) cells
- Molecular mass:  $\approx$  22 kDa (SDS-PAGE)
- Chromosome: 9
- UniProt ID: P01566
- Alternate name: IFN-alpha C
- Formulation: Recombinant human IFN- $\alpha$ 10 was lyophilized from a 0.2  $\mu$ m filtered phosphate buffer solution (pH 7.4) containing 2% human serum albumin (HSA) and 5% saccharose.
- Solubility: 100  $\mu$ g/ml in water

## APPLICATIONS

- Cellular assays (e.g. using HEK-Blue™ IFN- $\alpha$ / $\beta$  cells)
- ELISA.

## PRODUCT DESCRIPTION

Natural human IFN- $\alpha$ 10 is induced in peripheral blood mononuclear cells upon incubation with CpG-oligonucleotides and in plasmacytoid dendritic cells upon incubation with CpG-oligonucleotides or imiquimod<sup>1</sup>.

This subtype of IFN- $\alpha$  is classified among the top three strongest ISG inducers<sup>2,3</sup> with high anti-viral activity *in vitro* against human metapneumovirus<sup>4</sup> and hepatitis C virus (HCV)<sup>5</sup>. Human IFN- $\alpha$ 10 displays a strong capacity to induce IFIT1, CXCL10, CXCL11, ISG15 and CCL8<sup>3</sup>. Human IFN- $\alpha$ 10 expression is IFN- $\alpha$  receptor-dependent as it is induced by other IFN- $\alpha$  subtypes upon infection with low doses of Sendai virus *in vitro*<sup>6</sup>.

InvivoGen provides a glycosylated recombinant human IFN- $\alpha$ 10 produced in CHO cells. Of note, glycosylation stabilizes proteins against physicochemical instabilities.

1. Hillyer P. *et al.*, 2012. Expression profiles of human interferon-alpha and interferon-lambda subtypes are ligand- and cell-dependent. *Immunol. Cell. Biol.* 90(8):774. 2. Kurunganti S. *et al.*, 2014. Production and characterization of thirteen human type-I interferon- $\alpha$  subtypes. *Protein Expr. Purif.* 103: 75. 3. Moll H.P. *et al.*, 2011. The differential activity of interferon- $\alpha$  subtypes is consistent among distinct target genes and cell types. *Cytokines.* 53:52. 4. Scagnolari C. *et al.*, 2011. In vitro sensitivity of human metapneumovirus to type I interferons. *Viral Immunol.* 24(2):159. 5. Koyama T. *et al.*, 2006. Divergent activities of interferon-alpha subtypes against intracellular hepatitis C virus replication. *Hepatology Res.* 34(1):41. 6. Zaritsky L.A. *et al.*, 2015. Virus multiplicity of infection affects type I interferon subtype induction profiles and interferon-stimulated genes. *J. Virol.* 89:11534.

## METHODS

### Preparation of stock solution (20 $\mu$ g/ml):

1. Add 50  $\mu$ l endotoxin-free water (provided) to 1  $\mu$ g of recombinant hIFN- $\alpha$ 10 and mix gently by pipetting until completely resuspended.
2. Use immediately or prepare aliquots and store at -80 °C for 4 months. Avoid freeze-thaw cycles.
3. Further dilutions can be prepared in the appropriate aqueous buffer, such as cell culture medium containing serum.

## RELATED PRODUCTS

Product	Description	Cat. Code
Anti-hIFN- $\alpha$ -IgG	Neutralizing antibody	mabg-hifna-3
HEK-Blue™ IFN- $\alpha$ / $\beta$ cells	Type I IFN reporter cells	hkb-ifnab
Quanti-Blue™ Detection	Detection reagent	rep-qbs
LumiKine™ Xpress hIFN- $\alpha$ 2.0	Bioluminescent ELISA kit	luex-hifnav2

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)