

Recombinant human IFN- α 1

Mammalian cell-expressed human interferon alpha 1 (alpha D) with HSA

Cat. code: rcyc-hifna1

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

For research use only

Version 23A16-NJ

PRODUCT INFORMATION

Contents

- 1 μ g of lyophilized recombinant human IFN- α 1
- Note: This product is sterile filtered prior to lyophilization.
- 1.5 ml endotoxin-free water

Storage and stability

- Recombinant human IFN- α 1 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- α 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/ μ g
- The biological activity has been confirmed using HEK-Blue™ IFN- α / β 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 21 kDa (SDS-PAGE)
- Chromosome: 9
- UniProt ID: P01562
- Alternate name: IFN-alpha D
- Formulation: Recombinant human IFN- α 1 was lyophilized from a 0.2 μ m filtered phosphate buffer solution (pH 7.4) containing 2% human serum albumin (HSA) and 5% saccharose.
- Solubility: 100 μ g/ml in water

APPLICATIONS

- Cellular assays (e.g. using HEK-Blue™ IFN- α / β cells)
- ELISA.

PRODUCT DESCRIPTION

Despite being encoded by distinct genes, human IFN- α 1 and human IFN- α 13 display identical protein sequences¹. Human IFN- α 1 is induced in peripheral blood mononuclear cells (PBMCs) upon incubation with CpG-oligonucleotides and in plasmacytoid dendritic cells upon incubation with CpG-oligonucleotides or imiquimod². Viral infection of PBMCs with herpes simplex virus (HSV), Newcastle disease virus (NDV), and respiratory syncytial virus (RSV) induce high production of IFN- α 1³.

This subtype of IFN- α is classified as the weakest ISG inducer^{1,4} with the lowest anti-viral activity in vitro against influenza A virus¹ and hepatitis C virus (HCV)⁵. Similar to IFN- α 2 and α -8 subtypes, human IFN- α 1 expression occurs early post-infection with Sendai virus in vitro, regardless of the viral multiplicity of infection (MOI) and may deliver a 'priming' signal to neighboring cells⁶.

InvivoGen provides a glycosylated recombinant human IFN- α 1 produced in CHO cells. Of note, glycosylation stabilizes proteins against physicochemical instabilities.

1. Moll H.P. et al., 2011. The differential activity of interferon- α subtypes is consistent among distinct target genes and cell types. *Cytokines*. 53:52. 2. 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. 3. Löseke S. et al., 2003. Differential expression of IFN- α subtypes in human PBMC: evaluation of novel real-time PCR assays. *J. Immunol. Methods*. 276(1-2):207. 4. Kurunganti S. et al., 2014. Production and characterization of thirteen human type-I interferon- α subtypes. *Protein Expr. Purif.* 103: 75. 5. George J. & Mattapallil J.J., 2018. Interferon- α subtypes as an adjunct therapeutic approach for human immunodeficiency virus functional cure. *Front. Immunol.* 9:999. 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 μ g/ml):

1. Add 50 μ l endotoxin-free water (provided) to 1 μ g of recombinant hIFN- α 1 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- α -IgG	Neutralizing antibody	mabg-hifna-3
HEK-Blue™ IFN- α / β cells	Type I IFN reporter cells	hkb-ifnab
Quanti-Blue™ Detection	Detection reagent	rep-qbs
LumiKine™ Xpress hIFN- α 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