

IFN α

Interferon-Alpha 1 human human, recombinant, *E. coli*

Cat. No.	Amount
PR-450	10 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in PBS pH 7.4.
For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

Molecular Weight

20 kDa

Purity

$\geq 95\%$ by SDS-PAGE and RP-HPLC

Description

Interferons are cytokines that are widely known to induce potent anti-viral activity. Interferon α exerts a variety of other biological effects, including antitumor and immunomodulatory activities and are increasingly used clinically to treat a range of malignancies, myelodysplasias and autoimmune diseases.

IFNA1 Human Recombinant produced in E.Coli is a single, non-glycosylated, polypeptide chain containing 167 amino acids and having a molecular mass of 19.5 kDa. The IFN-alpha 1 is purified by proprietary chromatographic techniques.

Selected References:

- Birk *et al.* (2006) IFN-alpha induces apoptosis of adipose tissue cells. *Biochem Biophys Res Commun.* **345**:669.
Luft *et al.* (2002) IFN-alpha enhances CD40 ligand-mediated activation of immature monocyte-derived dendritic cells. *Int Immunol.* **14**:367.
Liu *et al.* (2001) High-yield expression and purification of human interferon alpha-1 in *Pichia pastoris*. *Protein Expr Purif.* **22**:381.