Recombinant Human IFN-α2b Protein Data Sheet

Catalog #	hRP-A0813-EF001
Size	25 μg
Protein Name	Human interferon alpha 2
Protein Symbol	IFN-α2b
Original Source	Homo sapiens
Expression System	E.coli
GenBank Accession #	NM 000605.2
Uniprot Accession #	P01563
Description	Interferon belongs to the cytokines family of proteins which are produced by mammalian cells in response to foreign agents such as viruses and tumor cells. It stimulates the production of two enzymes protein kinase and an oligoadenylate synthetase. Interferon alpha is a type 1 class of interferon produced by human leukocytes and has antiviral properties. Interferon alpha 2b is a recombinant protein produced in E. coli with a molecular mass of 19269 Dalton. It is a single, non-glycosylated, polypeptide chain containing 165 amino acids. The Interferon-alpha 2b gene was derived from human leukocytes. INF alpha 2b protein was purified using affinity chromatography.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	nature
Peptide Length	188aa
Molecular Weight	21.6kDa
pI	6.8
Activity	NA
Storage	Storage buffer: PBS containing 0.1% BSA.Store at -80°C and avoid repeated freeze-thaw cycles.
Reference:	 Immunogenicity of interferon-alpha 2 in therapy: structural and physiological aspects Structural and biological characterization of pegylated recombinant interferonalpha-2b and its therapeutic implications Zinc mediated dimer of human interferon-alpha 2b revealed by X-ray crystallography Effect of alpha-interferon treatment in patients with hepatitis B e antigen-positive chronic hepatitis B. A meta-analysis interferon alpha-2a (40 kDa) in the treatment of chronic hepatitis B. Interferon-based treatment of chronic hepatitis C



GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101 Rockville, MD 20850

USA

Phone: 301-762-0888 Toll free: 1-301-762-0888 Fax: 301-762-3888

Web: www.genecopoeia.com
Inquiry: inquiry@genecopoeia.com

Technical Support: support@genecopoeia.com