

# PRODUCT INFORMATION



## IFN- $\beta$ (human, recombinant)

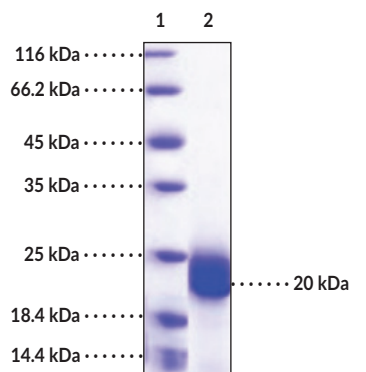
Item No. 31528

### Overview and Properties

**Synonyms:** Fibroblast Interferon, IFNB1, Interferon- $\beta$   
**Source:** Active recombinant human IFN- $\beta$  expressed in CHO cells  
**Amino Acids:** 22-187  
**Uniprot No.:** P01574  
**Molecular Weight:** 20 kDa  
**Storage:** -80°C (as supplied)  
**Stability:**  $\geq 1$  year  
**Purity:**  $\geq 95\%$  estimated by SDS-PAGE  
**Supplied in:** Lyophilized from sterile 10 mM sodium acetate with 40 mM Arg, 0.12 mg/ml Met, 146 mg/ml sucrose, and 0.1% PF-68, pH 4.5  
**Endotoxin Testing:**  $< 1.0$  EU/ $\mu$ g, determined by the LAL endotoxin assay  
**Bioactivity:** See figures for details

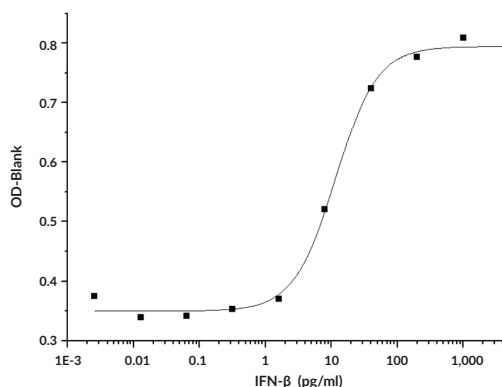
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: IFN- $\beta$

SDS-PAGE Analysis of IFN- $\beta$ .



IFN- $\beta$  Bioactivity in an Antiviral Assay Using WISH Cells Infected with Vesicular Stomatitis Virus. The EC<sub>50</sub> value for this effect is typically 2-20 pg/ml.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/23/2020

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Interferon- $\beta$  (IFN- $\beta$ ) is a cytokine and type I IFN with roles in antiviral responses and regulation of innate and adaptive immunity.<sup>1</sup> It is produced mainly by fibroblasts in response to viral pathogens, which are detected by a diverse repertoire of pathogen recognition receptors (PRRs).<sup>1,2</sup> IFN- $\beta$  binds to the IFN- $\alpha/\beta$  receptor (IFNAR) and induces signal transduction through the canonical JAK/STAT signaling pathway to induce the expression of IFN-stimulated genes (ISGs), which encode proteins that have antiviral, antiproliferative, or immunomodulatory properties, leading to the induction of an antiviral state in infected and neighboring cells and inhibiting viral replication.<sup>1</sup> IFN- $\beta$  inhibits the replication of a variety of viral pathogens, including severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2), the causative agent of COVID-19, and induces cell cycle arrest at the S phase and apoptosis in Huh7 hepatocellular carcinoma cells *in vitro*.<sup>3,4</sup> Formulations containing IFN- $\beta$  have been used in the treatment of multiple sclerosis. Cayman's IFN- $\beta$  (human, recombinant) protein can be used for cell-based assays. This protein consists of 166 amino acids, has a calculated molecular weight of 20 kDa, and a predicted N-terminus of Met22 after signal peptide cleavage.

## References

---

1. Ivashkiv, L.B. and Donlin, L.T. Regulation of type I interferon responses. *Nat. Rev. Immunol.* **14**(1), 36-49 (2014).
2. Haller, O., Kochs, G., and Weber, F. The interferon response circuit: Induction and suppression by pathogenic viruses. *Virology* **344**(1), 119-130 (2006).
3. Mantlo, E., Bukreyeva, N., Maruyama, J., *et al.* Antiviral activities of type I interferons to SARS-CoV-2 infection. *Antiviral Res.* **179**, 104811 (2020).
4. Murata, M., Nabeshima, S., Kikuchi, K., *et al.* A comparison of the antitumor effects of interferon- $\alpha$  and  $\beta$  on human hepatocellular carcinoma cell lines. *Cytokine* **33**(3), 121-128 (2006).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM