

IFN- α -2b (human), (recombinant) (active)

Active enzyme with antiviral and antiproliferative properties.

IFN- α -2 is a cytokine that displays antiviral and antiproliferative properties. It is comprised of three major subtypes, which include IFN- α -2a, IFN- α -2b, and IFN- α -2c. IFN- α -2b is the predominant subtype in the human genome. It binds to its receptor and activates the JAK-STAT pathway, which stimulates expression of the host genes responsible for antiviral activity. IFN- α -2b is used to treat many diseases, such as hepatitis B and C, melanoma, Kaposi's sarcoma, chronic myeloid lymphoma, and angioblastoma. The combination of IFN- α -2b with ribavirin, a viral inhibitor, has become the standard treatment for chronic hepatitis C virus infection.

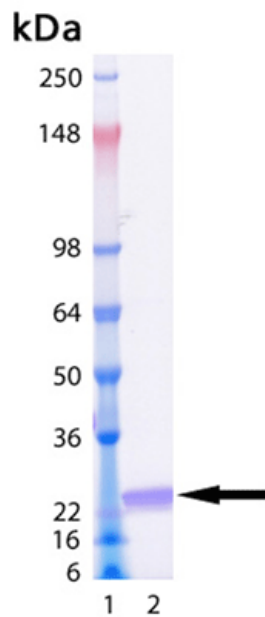
Ordering Information

[Order Online »](#)

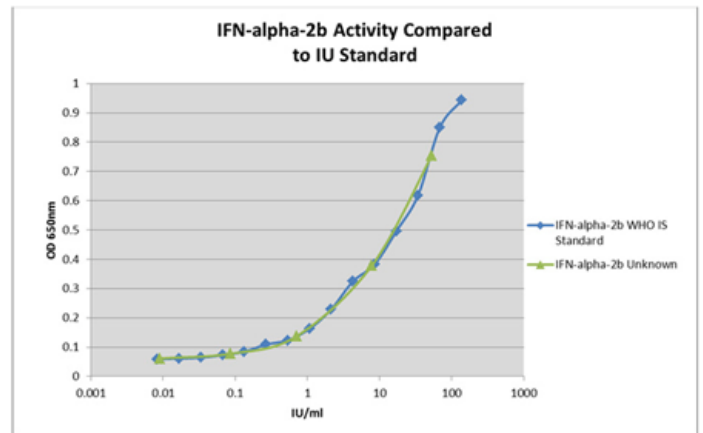
ENZ-PRT192-0010	10 μ g
ENZ-PRT192-5010	5x10 μ g

Manuals, SDS & CofA

[View Online »](#)



Coomassie stained SDS-PAGE. Lane 1, Molecular weight marker. Lane 2, 1.0µg Human IFN- α -2b (ENZ-PRT192).



ENZ-PRT192 activity was determined by comparing against the WHO International Standard for Interferon Alpha 2b (NIBSC, Code 95/566) using the HEK-Blue™ hTLR5 reporter cell line (Invivogen). Briefly, cells were plated at 50,000 cells per well in 100µl of test media. Known amounts of WHO standard and serially diluted ENZ-PRT192 were added in a volume of 100µl. The cells were grown for 20-24 hours at 37°C + 5% CO₂. After incubation, 180µl of Quanti-Blue™ was added to 20µl of HEK-Blue™ hTLR5 reporter cell supernatant and the color was allowed to develop for 3 hours at 37°C + 5% CO₂. Activity was determined by interpolating the OD at 650nm of the ENZ-PRT192 titer off of the WHO standard curve.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	>5000 IU/μg
Alternative Name	Interferon-α-2b
Concentration	0.1 mg/ml when reconstituted as suggested.
Endotoxin Content	<0.1EU/μg protein (LAL test)
Formulation	Lyophilized. In PBS.
MW	~21kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Affinity purified.
Reconstitution	Reconstitute with 100μl sterile distilled water.
Source	Produced in HEK cells. The human cytokine (aa 24-188) is fused at the N-terminus to a FLAG®-tag followed by an 8 AA linker.
Technical Info / Product Notes	FLAG® is a registered trademark of Sigma-Aldrich Co.
UniProt ID	P01563



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com