

Human Interferon Beta 1a, mammalian

Citations: 5

[View Online »](#)

Ordering Information

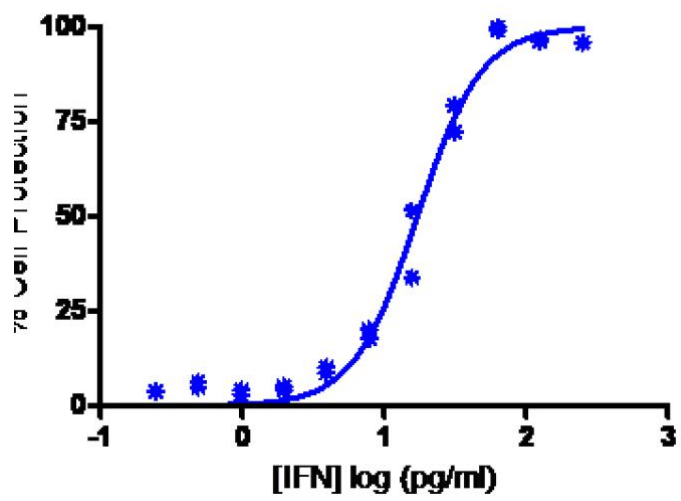
[Order Online »](#)

PBL-11415-1	0.1MU
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Human IFN Beta 1a, mammalian, recombinant protein expressed in CHO cells
- Supplied frozen with a carrier protein for retention of full activity
- Used for MS research studies



Titration of Human IFN-Beta in A549 Cells

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon-β1a, IFN-β1a

Endotoxin Content < 1 EU/μg

Formulation Liquid. In 50mM sodium acetate, pH 5.5, containing 0.1% BSA.

MW ~23kDa

Purity ≥95%

Purity Detail Proprietary chromatography methods

Quantity 0.1ml

Source Produced in CHO cells.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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