

# Human Interferon Beta 1a, mammalian

Citations: 4

[View Online »](#)

## Ordering Information

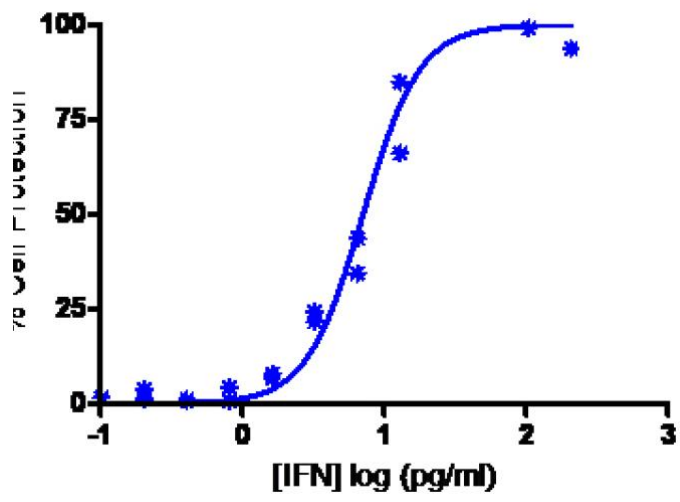
[Order Online »](#)

PBL-11410-2	1MU
-------------	-----

## Manuals, SDS & CofA

[View Online »](#)

- Human IFN Beta 1a, mammalian, recombinant protein expressed in CHO cells
- Supplied without a carrier protein and as a liquid to maintain higher specific activity
- Used for MS research studies



Titration of mammalian expressed Human IFN-Beta on A549 Cell Line

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Interferon- $\beta$ 1a, IFN- $\beta$ 1a
Endotoxin Content	< 1 EU/ $\mu$ g
Formulation	Liquid. In 10mM acetic acid.
MW	~23kDa
Purity	$\geq$ 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)