

# Mouse Interferon Beta

Citations: 4

[View Online »](#)

## Ordering Information

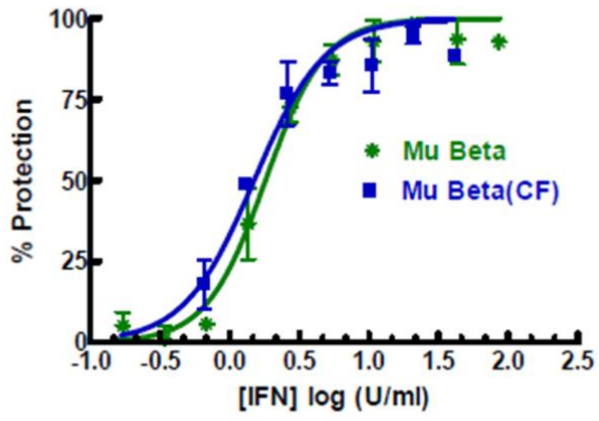
[Order Online »](#)

PBL-12400-1	0.1MU
-------------	-------

## Manuals, SDS & CofA

[View Online »](#)

- E.coli expressed Mouse IFN Beta protein with a single non-glycosylated polypeptide chain
- Increased stability with engineered disulfide bond
- 95% or greater purity



Antiviral Activities of Mouse IFN-Beta (112400) and carrier-free Mouse IFN-Beta (12401) on Mouse L929 Cells infected with EMCV

## Handling & Storage

<b>Use/Stability</b>	Thaw product vial by incubation in cold tap water until just thawed. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA).
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Interferon- $\beta$ , IFN- $\beta$
<b>Endotoxin Content</b>	< 1 EU/ $\mu$ g
<b>Formulation</b>	Liquid. In PBS containing 0.1% BSA.
<b>MW</b>	~19.7kDa
<b>Purity</b>	$\geq$ 95%
<b>Purity Detail</b>	A combination of ion exchange, hydrophobic interaction and size exclusion chromatography.
<b>Quantity</b>	0.1ml
<b>Source</b>	Produced in <i>E. coli</i> .



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)