

# Anti-Mouse Interferon Beta, Rabbit IgG, Protein A purified (PAb)

Binds to and neutralizes mouse IFN beta

## Ordering Information

[Order Online »](#)

PBL-32401-1	2x10kU
-------------	--------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	IFN- $\beta$ , Interferon- $\beta$
Formulation	Liquid. In 0.3 M NaHCO <sub>3</sub> ; 0.2 M NaCl
Host	Rabbit
Immunogen	Mouse IFN- $\beta$ (interferon- $\beta$ ).
Species Reactivity	Mouse



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)