

# Anti-Human IFN Beta Antibody, Goat IgG, Affinity Purified, neutralizing (PAb)

Neutralizes human interferon beta

## Ordering Information

[Order Online »](#)

PBL-31420-1	100µg
-------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	IFN- $\beta$ , Interferon- $\beta$
Application	ELISA, WB
Crossreactivity	Does not cross-react with IFN- $\alpha$ or IFN- $\gamma$ .
Formulation	Lyophilized from a 0.2 $\mu$ m filtered solution in phosphate – buffered saline (PBS) with 5% trehalose
Host	Goat
Immunogen	Recombinant human IFN- $\beta$ .
Species Reactivity	Human



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)