

# Anti-Mouse Interferon Gamma, Clone RMMG- 1 (MAb)

Neutralizes mouse interferon gamma; binds to mouse interferon gamma with high affinity

## Ordering Information

[Order Online »](#)

PBL-22500-1	500µg
-------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	IFN- $\gamma$ , Interferon- $\gamma$
Application	ELISA, WB
Clone	RMMG-1
Formulation	Liquid. In PBS containing 0.1% BSA.
Host	Rat
Immunogen	Mouse IFN- $\gamma$ (interferon- $\gamma$ ).
Isotype	IgG1
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)