

# Anti-Human Interferon Gamma, Clone MMHG- 2 (MAb)

Binds to but does not neutralize human interferon gamma

## Ordering Information

[Order Online »](#)

PBL-21550-1	100µg
-------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	IFN- $\gamma$ , Interferon- $\gamma$
<b>Clone</b>	MMHG-2
<b>Formulation</b>	Liquid. In PBS containing 0.1% BSA.
<b>Host</b>	Mouse
<b>Immunogen</b>	Human IFN- $\gamma$ (interferon- $\gamma$ ).
<b>Isotype</b>	IgG1
<b>Species Reactivity</b>	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)