

Anti-Human Interferon Gamma, Clone MMHG- 1 (MAb)

Binds to and neutralizes human interferon gamma

Ordering Information

[Order Online »](#)

PBL-21500-1	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- γ , Interferon- γ
Clone	MMHG-1
Formulation	Liquid. In PBS containing 0.1% BSA.
Host	Mouse
Immunogen	Human IFN- γ (interferon- γ).
Isotype	IgG1
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)