

Anti-Mouse Interferon Alpha, Chicken IgY, Purified (PAb)

Binds to but not neutralize mouse IFN alpha

Ordering Information

[Order Online »](#)

PBL-32120-1	500µg
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	After reconstitution, 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- α , Interferon- α
Application	WB
Formulation	Liquid. In PBS containing 0.1% BSA.
Host	Chicken
Immunogen	Recombinant mouse IFN- α (interferon- α).
Species Reactivity	Mouse



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com