

Anti-Human Interferon Alpha, Rabbit Serum (PAb)

Binds well to human IFN alpha A (2a), A/D and K (6)

Ordering Information

[Order Online »](#)

PBL-31101-1	0.1MU
-------------	-------

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- α , Interferon- α
Formulation	Neat serum in PBS.
Host	Rabbit
Immunogen	Recombinant human IFN- α A (interferon- α A).
Species Reactivity	Human



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