

Anti-Human Interferon Alpha (All Subtypes), Rabbit Serum (PAb)

Binds to and blocks biological activity of all human interferon alpha subtypes

Ordering Information

[Order Online »](#)

PBL-31135-1	1MU
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

- Rabbit polyclonal antibody that neutralizes all Human IFN Alpha subtypes (IFN-alpha 2a, IFN-alpha 2b, IFN-alpha 8, IFN-alpha 10, IFN-alpha 1, IFN-alpha 21, IFN-alpha 5, IFN-alpha 14, IFN-alpha 17, IFN-alpha 7, IFN-alpha 6, IFN-alpha D, IFN-alpha M1, IFN-alpha 4, IFN-alpha 16)
- Anti-human polyclonal antibody designed to accurately bind with and block biological activity of all human interferon alpha subtypes

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN alpha

Formulation Neat serum diluted in phosphate buffered saline (PBS)

Host Rabbit

Immunogen Mixture of all recombinant human interferon alpha subtypes (IFNA2, IFNA8, IFNA10, IFNA1, IFNA21, IFNA5, IFNA14, IFNA17, IFNA7, IFNA6, IFNA4, and IFNA16)

Species Reactivity Human

Specificity Binds to and blocks biological activity of all human interferon alpha subtypes



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