

Anti-Human Interferon Alpha, Sheep Serum (PAb)

Ordering Information

[Order Online »](#)

PBL-31110-1	0.1MU
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Neutralizes all Human IFN Alpha subtypes
- Suitable for Direct Binding ELISA

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 Sheep

Immunogen Mixture of recombinant human IFN 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