

Human Type I Interferon Neutralizing Antibody Mixture

Mixture of monoclonal and polyclonal antibodies directed against human Type I interferon receptor and several Type I interferons

Citations: 5

[View Online »](#)

Ordering Information

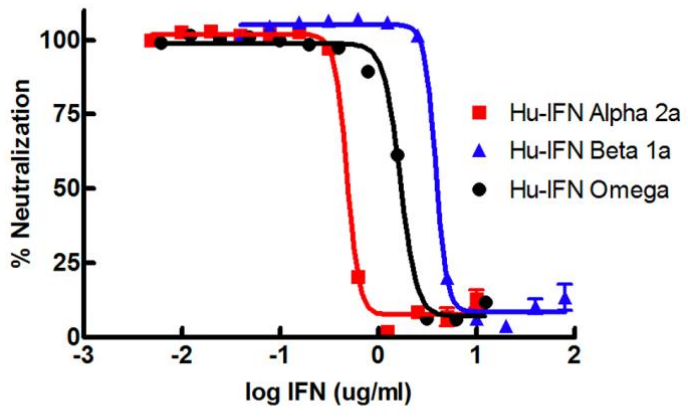
[Order Online »](#)

PBL-39000-1	0.2ml
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Mixture of monoclonal and polyclonal antibodies directed against human Type 1 Interferons (IFNs) and Type 1 Receptor Subunit 2 (IFNAR2)
- Effectively **neutralizes** biological activity of human IFN-alpha, IFN-beta, IFN-omega, IFN-kappa, and IFN-epsilon
- Improves **neutralization** of multiple Type I IFNs often found in complex samples by employing a broadly neutralizing custom antibody cocktail
- Unconjugated



Human Type I IFN Neutralization in A549 Cells

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN

Crossreactivity No cross reactivity observed with the following human interferons: gamma, lambda 1, lambda 2, lambda 3; or mouse interferons: alpha, beta, gamma

Formulation Supplied frozen in phosphate-buffered saline (PBS)

Purity Detail Purified by a combination of Protein G affinity chromatography and other conventional chromatographical methods.

Species Reactivity Human

Specificity Blocks biological activity of human Type I interferons (alpha, beta, omega, kappa, epsilon)



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