

Anti-Human Interferon Alpha/Beta Receptor Chain 2, Clone MMHAR-2 (MAb)

Neutralizes human interferon alpha receptor; interacts with extracellular domain; binds to human interferon alpha receptor with high affinity; blocks biological action of Type I IFNs

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

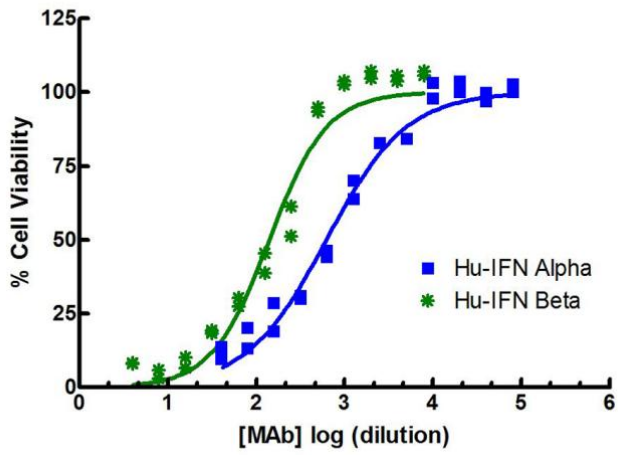
PBL-21385-1	50µg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

- Anti-IFNAR2 antibody (clone MMHAR-2) to human Interferon (Alpha, Beta and Omega) Receptor 2
- IFN Receptor antibody blocks the biological action of Human Type I Interferons
- Suitable for use in Type I IFN neutralization, flow cytometry, immunohistochemistry, and immunoprecipitation

**Mouse MAb to Human
IFN Alpha/Beta Receptor 2 (MMHAR-2)**



Neutralization Effect of PBL Anti-Human IFNAR2
Monoclonal Antibody (21385) on Human IFN-Alpha 2a
and IFN-Beta Activities in an A549 Cell/EMCV CPE
assay

Handling & Storage

Handling	For long term storage keep unconjugated antibody at -20°C. Avoid freeze/thaw cycles. Keep conjugated formats at +4°C. Do not freeze. Protect from light.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	IFN- α/β Receptor Chain 2, Interferon- α/β Receptor Chain 2
Clone	MMHAR-2
Formulation	Liquid. PBL-21385-1 (Unlabelled): In PBS containing 0.1% BSA. PBL-21385-3 (R-PE labelled): In PBS, pH 7.4, containing 1.5% BSA and 0.1% Kathon as preservative.
Host	Mouse
Immunogen	Human IFN alpha/beta RC2
Isotype	IgG2a
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