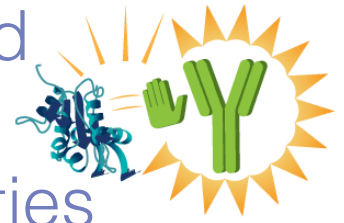




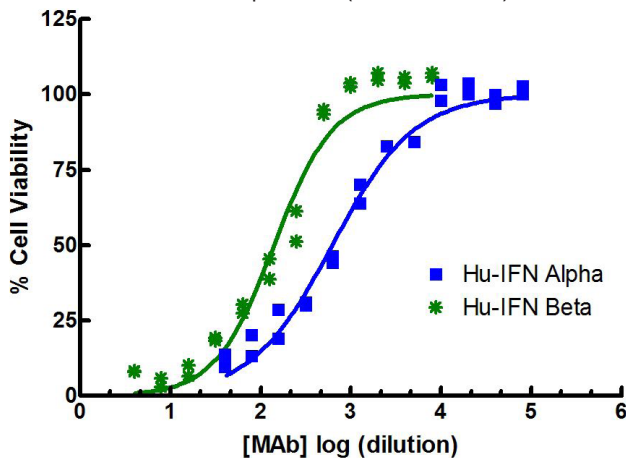
Select PBL antibodies and proteins to neutralize human & mouse IFN activities



21385-1

Anti-Human Interferon Alpha/Beta Receptor Chain 2 (IFNAR2), Clone MMHAR-2 (MAb)

Figure 1. Mouse MAb to Human IFN α/β Receptor 2 (MMHAR-2)



- Neutralizes biological activity of Human Type I IFNs (α , β , and ω)
- Blocks Type I interferon receptor complex
- Suitable for use in Type I IFN neutralization, flow cytometry, immunohistochemistry, and immunoprecipitation

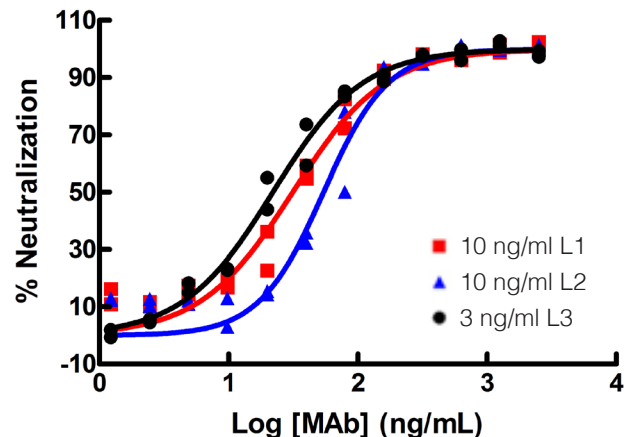
21885-1

Anti-Human Interferon Lambda Receptor 1, Clone MMHLR-1 (MAb)

- Neutralizes human interferon lambda receptor 1

Sample	EC50 (ng/ml)
10 ng/ml L1	31.3
10 ng/ml L2	54.3
3 ng/ml L3	22.0

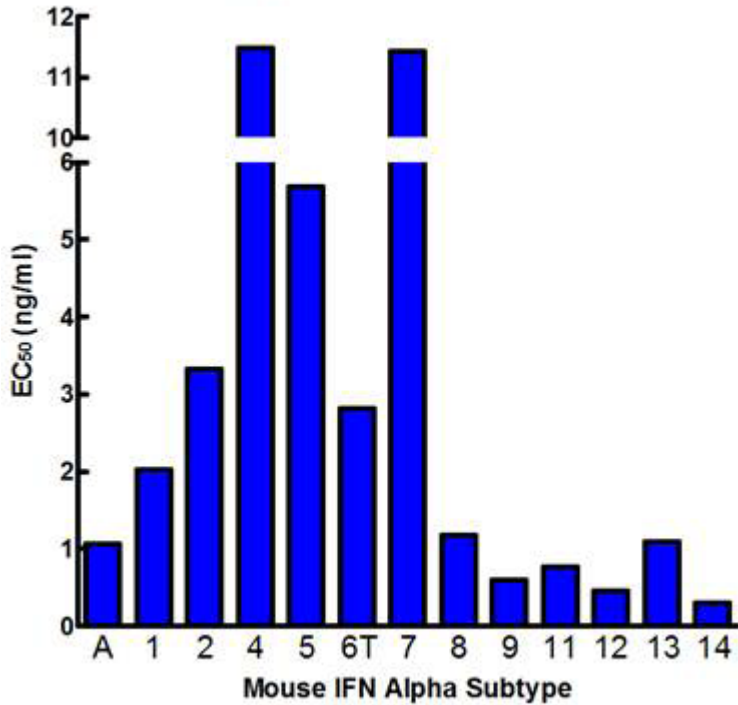
Figure 2. Neutralizing MAb against IFNLR1



12185-1

C12R Mouse Interferon Alpha Decoy Receptor

Figure 3. C12R Neutralization of Mouse IFN α Subtypes on L929 Cells



- Viral protein C12R neutralizes all 14 mouse IFN α subtypes *in vitro*
- Neutralizes all human IFN α subtypes

31110-1

Anti-Human Interferon Alpha, Sheep Serum (PAb)

Figure 4. Neutralization of Human IFN α subtypes on A549 cells

- Neutralizes all human interferon alpha subtypes
- Suitable for use in neutralization and Direct Binding ELISA

