



Quality and Value - A Perfect Dimer!

High quality Interferon Gamma products from PBL Assay Science:

- High Specific Activity
- Low in aggregates
- > 98% dimerized

Did you know that Interferon Gamma is only active as a protein dimer?

Interferon Gamma are members of the Type II class of interferons. Type II IFN in humans is limited to the single Interferon Gamma gene. This IFN binds to the Type II receptor comprised of IFNGR1 and IFNGR2 chains. Interferon Gamma is produced by cells of the immune system such as B-cells, T-cells and NK cells. Interferon Gamma is glycosylated in mammalian cells and the active form functions as a homodimer. All of PBL's Interferon Gamma proteins are guaranteed dimers for maximum activity.

Figure 1. Chromatographic profile of IFN- γ showing dimerization and low aggregation

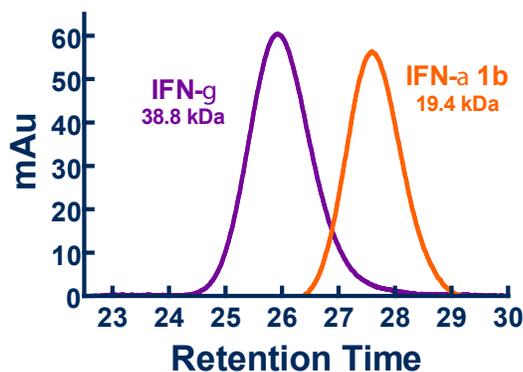
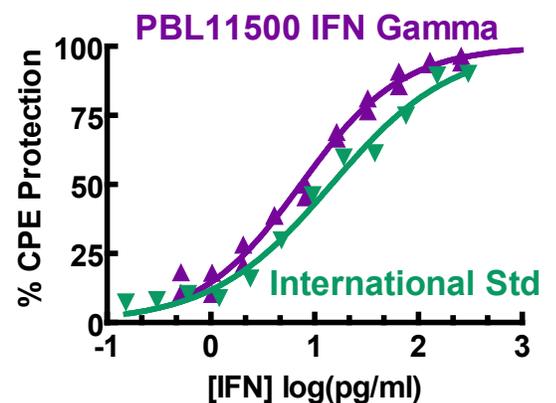


Figure 2. PBL IFN- γ and international gamma standard exhibit comparable antiviral activities



Product Information

Interferon Gamma Products		
11500-1	Human IFN Gamma	50 µg
11500-2	Human IFN Gamma	100 µg
13500-1	Rat IFN Gamma	10 µg
21500-1	Anti-Human IFN Gamma, Clone MMHG-1 (MAb)	500 µg
21550-1	Anti-Human IFN Gamma, Clone MMHG-2 (MAb)	100 µg
21585-1	Anti-Human IFN Gamma Receptor Chain 2, Clone MMHGR-2 (MAb)	250 µg
22500-1	Anti-Mouse IFN Gamma, Clone RMMG-1 (MAb)	500 µg
31500-1	Anti-Human IFN Gamma, Rabbit Serum (PAb)	2 x 10 ⁴ U
32500-1	Anti-Mouse IFN Gamma, Rabbit Serum (PAb)	2 x 10 ⁴ U
41500-1	VeriKine Human IFN Gamma ELISA Kit	1 plate
41580-1	NEW VeriKine Human IFN Gamma Receptor 1 ELISA Kit	1 plate

In addition to the standard products listed, PBL Assay Science offers customized production of cytokines, interferons, and antibodies. For details regarding these services please contact our customer service department by telephone (732-777-9123) or e-mail (info@pblassaysci.com).

