



Mouse Interferon Reagents and ELISAs

- High quality recombinant proteins for research
- Antibodies validated for biological assay use
- ELISAs for immunological assay development

As the drug discovery pathway shifts toward specifically targeted pathways and molecules, model systems continue to increase in importance. The ease of use and versatility of mouse models in the study of human cancer, infectious disease and autoimmunity provides understanding into the etiology of these diseases. Data obtained from studies using mouse models can provide insight into potential clinical side effects and the pathology of diseases.

PBL Assay Science provides a comprehensive mouse product line manufactured to the highest standards of quality, consistency and purity. PBL's products are designed to meet your scientific needs by providing superior quality and performance to ensure reliable and reproducible results.

Figure 1. Comparison of Mu Alpha A and Mu Alpha 4 Antiviral Activity

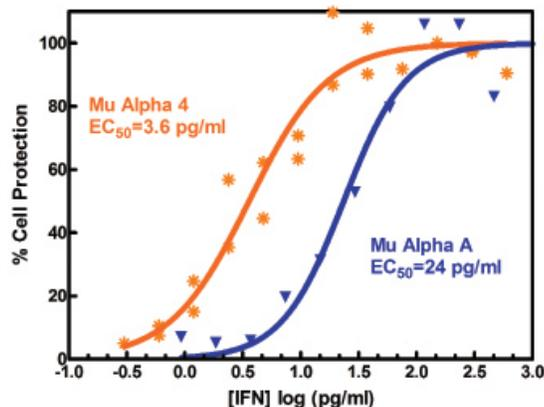
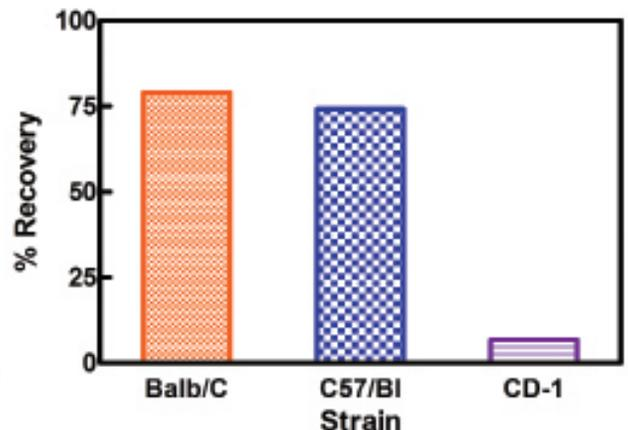


Figure 2. Recovery of Mu Alpha A in 50% Serum in Mu Alpha ELISA (42120)



Product Information

Interferons

12100-1	Mouse IFN Alpha A	1 x 10 ⁵ U
12105-1	Mouse IFN Alpha 1	1 x 10 ⁵ U
12115-1	Mouse IFN Alpha 4, mammalian	1 x 10 ⁵ U
12125-1	Mouse IFN Alpha 11	1 x 10 ⁵ U
12130-1	Mouse IFN Alpha 13	1 x 10 ⁵ U
12185-1	C12R Mouse IFN Alpha Decoy Receptor, neutralizing	10 µg
12185-2	C12R Mouse IFN Alpha Decoy Receptor, neutralizing	100 µg
12400-1	Mouse IFN Beta	1 x 10 ⁵ U
12401-1	Mouse IFN Beta (carrier-free)	1 x 10 ⁵ U
12405-1	Mouse IFN Beta, mammalian	1 x 10 ⁵ U
12410-1	Mouse IFN Beta, mammalian (carrier-free)	1 x 10 ⁶ U
12820-1	Mouse Interleukin-28b/IFN Lambda 3	25 µg
12821-1	Mouse Interleukin-28b/IFN Lambda 3 (carrier-free)	25 µg

- Recombinant proteins derived from bacterial or mammalian cells
- Consistent performance and biological activity
- > 95% purity
- Endotoxin level <1 EU/µg
- Suitable for use in a range of applications such as ELISAs and cytopathic effect assays
- Viral C12R IFN Antagonist neutralizes all mouse IFN alpha subtypes

Antibodies

22100-1	Anti-Mouse IFN Alpha, Clone RMMA-1 (MAb)	250 µg
22100-3	FITC Conjugated Anti-Mouse IFN Alpha, Clone RMMA-1 (MAb)	25 µg
22400-1	Anti-Mouse IFN Beta, Clone RMMB-1 (MAb)	250 µg
22500-1	Anti-Mouse IFN Gamma, Clone RMMG-1 (MAb)	500 µg
32100-1	Anti-Mouse IFN Alpha, Rabbit Serum (PAb)	2 x 10 ⁴ U
32120-1	Anti-Mouse IFN Alpha, Chicken IgY, Purified (PAb)	500 µg
32400-1	Anti-Mouse IFN Beta, Rabbit Serum (PAb)	2 x 10 ⁴ U
32401-1	Anti-Mouse IFN Beta, Rabbit IgG, Protein A Purified (PAb)	2 x 10 ⁴ U
32500-1	Anti-Mouse IFN Gamma, Rabbit Serum (PAb)	2 x 10 ⁴ U

- High specificity
- Monoclonal antibodies have > 95% purity
- Suitable for use in a range of applications such as ELISAs, neutralization, flow cytometry, and western blotting

ELISAs

42120-1	VeriKine Mouse IFN Alpha ELISA Kit	1 plate
42120-2	VeriKine Mouse IFN Alpha ELISA Kit	5 plate
42115-1	VeriKine-HS Mouse IFN Alpha All-Subtype ELISA Kit	1 plate
42400-1	VeriKine Mouse IFN Beta ELISA Kit	1 plate
42400-2	VeriKine Mouse IFN Beta ELISA Kit	5 plate
42410-1	VeriKine-HS Mouse IFN Beta Serum ELISA Kit	1 plate
42410-2	VeriKine-HS Mouse IFN Beta Serum ELISA Kit	5 plate
52500-1	VeriPlex Mouse Cytokine 9-Plex ELISA Kit	1 plate

- Mouse Alpha ELISAs provide global interferon alpha measurement in serum, plasma, and tissue culture media
- < 10% inter- and intra-assay CVs
- High sensitivity ELISA options provide accurate results ~2 hours

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).