



Mouse Interferon Products

Highly Active Mouse Interferon Proteins

High quality recombinant proteins derived from bacterial or mammalian cells

- Consistent performance and biological activity
- 95% or greater purity
- Endotoxin level < 1 EU/μg
- Suitable for use in a range of applications including ELISA & CPE assays

Featured:

Mouse IFN-Alpha 4, mammalian (Cat. No. 12115)

Recombinant Type I interferon exhibiting early induction

- Provides a better model of native mouse alpha 4 through authentic sequencing
- May play a role in preparing cells to express high levels of other mouse IFN-Alpha subtypes upon viral infection

C12R Mouse IFN-Alpha Decoy Receptor (Cat. No. 12185)

Viral C12R IFN antagonist selectively neutralizes all mouse IFN-Alpha subtypes

- No cross-reactivity with mouse IFN-Beta

Figure 1. (top right) Antiviral Activity Comparison of Mouse IFN-Alpha A and IFN-Alpha 4 (mammalian). Mouse IFN-Alpha 4 exhibits higher antiviral potency than Mouse IFN-Alpha A in this standard CPE assay.

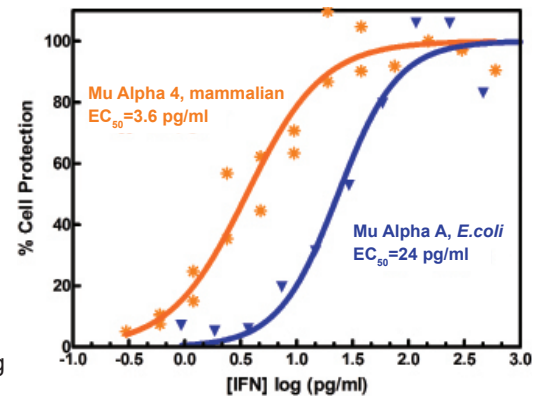
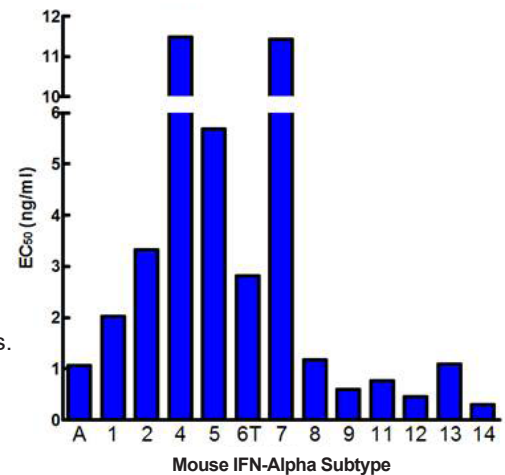


Figure 2. (bottom right) C12R Neutralization of Mouse IFN-Alpha Subtypes on L929 Cells.



Cat. No.	Description	Size
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, 2	C12R Mouse IFN-Alpha Decoy Receptor, neutralizing	10 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

To learn more, please contact PBL Assay Science or visit PBL's website.

Mouse Interferon Alpha (IFN- α) ELISAs

Achieve global detection and quantitation of all Mouse IFN-Alpha subtypes in sera, plasma, or Tissue Culture Media (TCM).

Featured:

Mouse IFN-Alpha All Subtype ELISA (Cat. No. 42115)

High Sensitivity

- Assay Standard Range: 2.38 – 152 pg/ml
- Recognizes all 14 IFN-Alpha subtypes
- Sensitive, low pg/ml global detection of Mouse IFN-Alpha subtypes
- Robust matrix tolerance including serum, plasma, and tissue culture media (TCM)
- Rapid ELISA provides total interferon alpha readout in ~ 2 hours
- High data reproducibility with < 10% inter- and < 8% intra-assay CVs

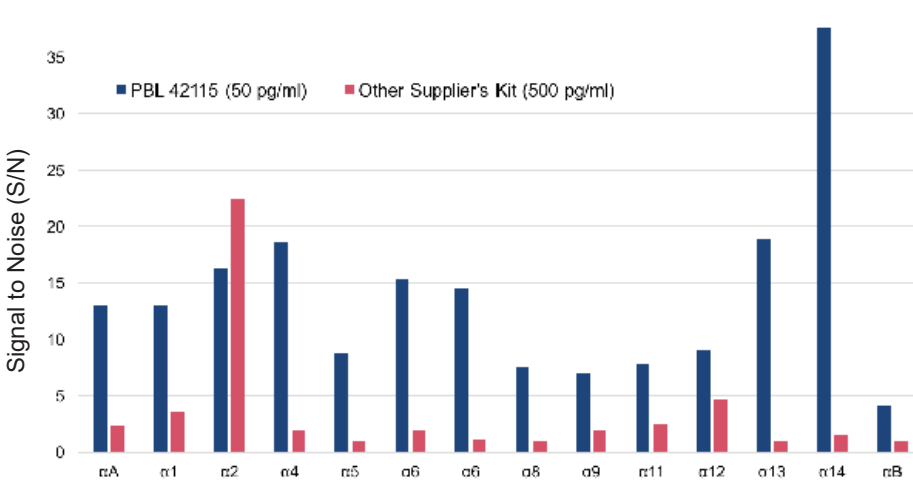


Figure 3. Signal to Noise Ratios of Mouse IFN-Alpha Subtypes in Mouse Serum.

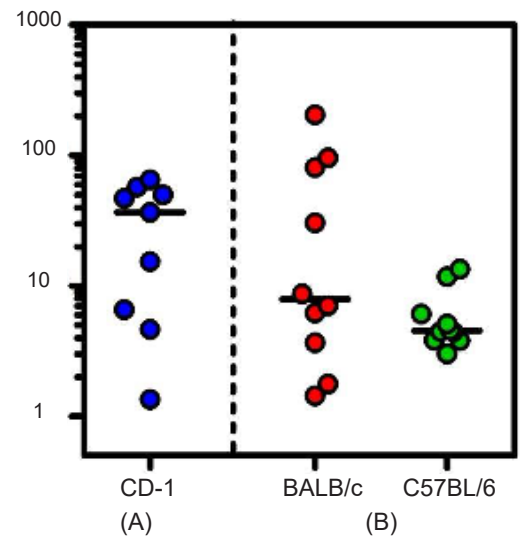


Figure 4. Levels of endogenous mouse IFN-Alpha quantified in pooled plasma (A) and individual sera (B).

	42120	42115
• Good •• Better ••• Best	Mu-IFN- α	Mu-IFN- α All Subtype
Global IFN-Alpha subtype measurement in serum or plasma	••	•••
Global IFN-Alpha subtype measurement in TCM	••	•••
Reproducible results with < 10% inter- & intra-assay CVs	•••	•••

Cat. No.	Description	Size
42120-1, 2	VeriKine Mouse IFN-Alpha ELISA Kit (TCM)	1 x 96-well plate, 5 x 96-well plate
42115-1	VeriKine-HS Mouse IFN-Alpha All Subtype ELISA Kit (S, P, TCM)	1 x 96-well plate

To learn more, please contact **PBL Assay Science** or visit **PBL's website**.

Mouse Interferon Beta (IFN-β) ELISAs

Superior detection and measurement of Mouse IFN-Beta in sera, plasma, or TCM at an unrivaled level of sensitivity.

Featured:

Mosue IFN-Beta Serum ELISA (Cat. No. 42410)

High Sensitivity

- Assay Standard Range: 0.94 – 60 pg/ml
- Accurately quantifies low levels of Mouse IFN-Beta in multiple matrices
- Provides > 90% spike-recovery in serum and plasma
- High precision assay reproducibility with ≤ 7% and ≤ 6% intra-assay CVs
- Total assay time of ~ 2 hours

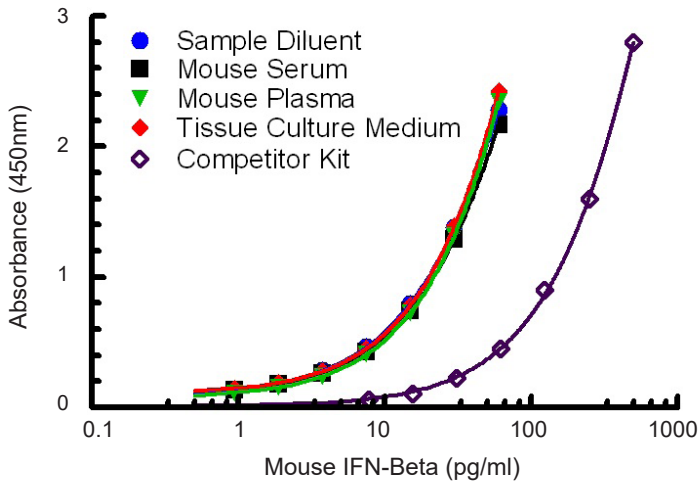


Figure 5. Representative Standard Curves. Competitor data obtained from published public source (provided for comparison use only).

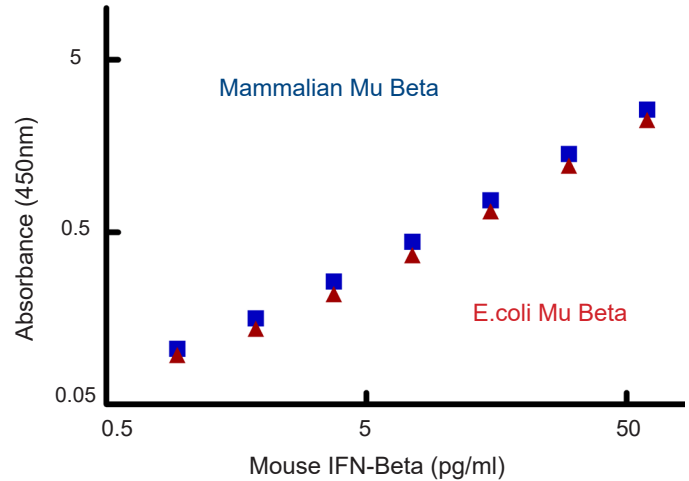


Figure 6. Comparison of ELISA Reactivity of Mammalian and E.coli-expressed Mouse IFN-Beta. Sensitive detection of mammalian and *E.coli*-expressed IFN-Beta.

	42400	42410
	Mu-IFN-β	Mu-IFN-β Serum
• Good •• Better ••• Best		
Accurate IFN-Beta measurement in serum or plasma		•••
Accurate IFN-Beta measurement in TCM	••	•••
Reproducible results with < 10% inter- & intra-assay CVs	••	•••

Cat. No.	Description	Size
42400-1, 2	VeriKine Mouse IFN-Beta ELISA Kit (TCM)	1 x 96-well plate, 5 x 96-well plate
42410-1, 2	VeriKine-HS Mouse IFN-Beta Serum ELISA Kit (S, P, TCM)	1 x 96-well plate, 5 x 96-well plate

To learn more, please contact PBL Assay Science or visit PBL's website.

Anti-Mouse IFN Monoclonal (MAb) & Polyclonal (PAb) Antibodies

Labeled & unlabeled antibodies for a variety of uses

- MAbs offer high specificity in applications where specific binding is desirable with > 95% purity
- PABs are preferable for neutralization because they target a variety of specific epitopes of a particular antigen

Featured:

Anti-Mouse IFN-Beta Antibody, Clone RMMB-1 (MAb) (Cat. No. 22400)

- Binds to mouse IFN-Beta (weakly neutralizes)
- Suitable for use in Direct Binding ELISA

Anti-Mouse IFN-Beta Antibody, Rabbit IgG, Protein A Purified (PAb) (Cat. No. 32401)

- Suitable for use in neutralization
- Suggested for use in ELISA, Western Blot, immunoprecipitation, and immunohistochemistry

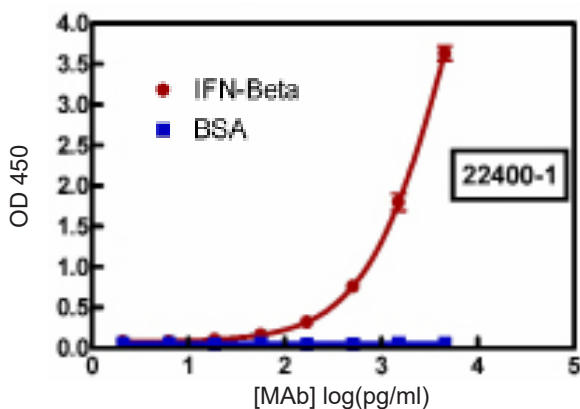


Figure 7. Direct Binding ELISA. Representative binding curve of Anti-Mouse IFN-Beta antibody to recombinant Mouse IFN-Beta and 1% BSA/ PBS.

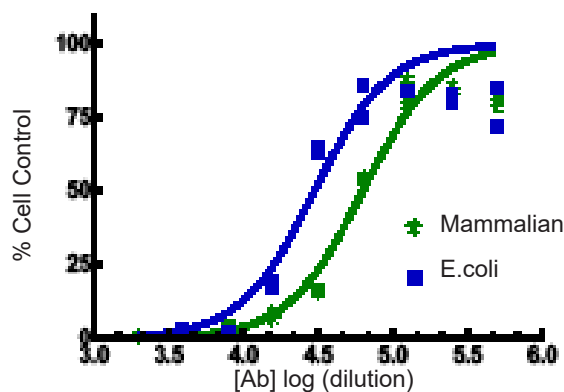


Figure 8. Titration against Mammalian IFN. Neutralization of 32401 Rabbit PAb stocks using mammalian-expressed Mouse IFN-Beta standard.

Cat. No.	Description	Size
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 units
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 units
32401-1	Anti-Mouse IFN-Beta, Rabbit IgG, Protein A Purified (PAb)	2 x 10 ⁴ units
32500-1	Anti-Mouse IFN-Gamma, Rabbit Serum (PAb)	2 x 10 ⁴ units

To learn more, please contact **PBL Assay Science** or visit **PBL's website**.



pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA
 T +1.732.777.9123 | F +1.732.777.9141 | E info@pblassaysci.com | W pblassaysci.com