

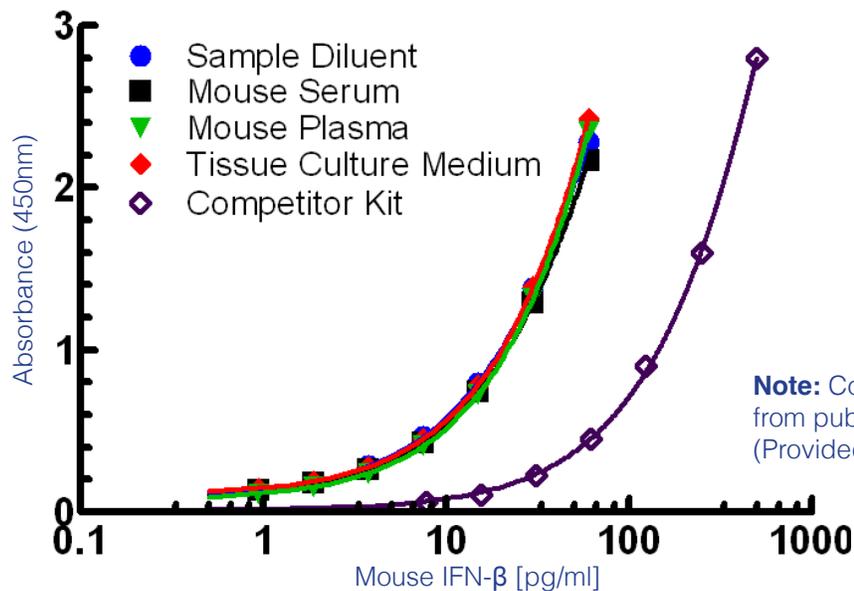


VeriKine-HS™ Mouse Interferon Beta Serum ELISA Kit

Catalog No. 42410

- The **world's most sensitive** Mouse IFN Beta ELISA
- Quantifies low level mouse IFN- β in **serum, plasma** and tissue culture media with an assay range of **0.94 - 60 pg/ml**
- Reliable serum and plasma spike recoveries of >90%
- Provides reproducible results with <10% CV
- Total assay time of **~ 2 hours**

Figure 1. Standard Curve



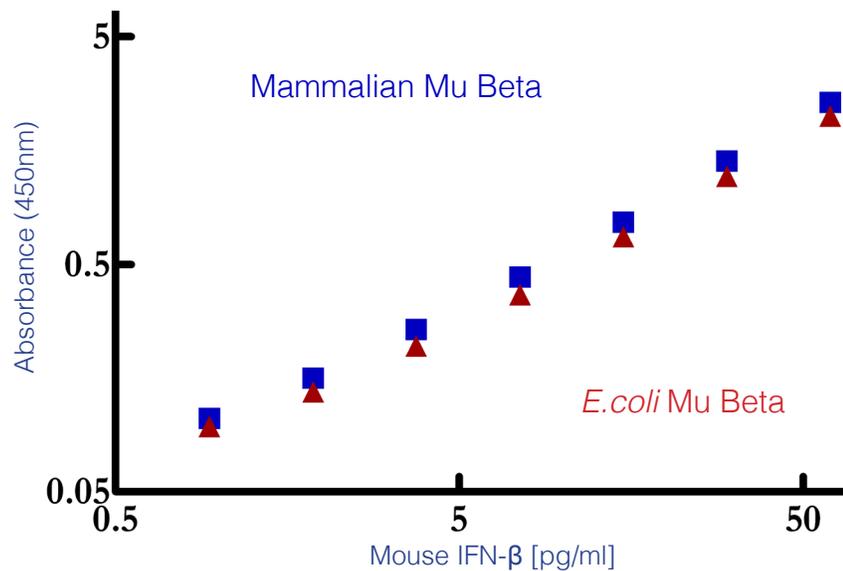
Note: Competitor data obtained from published public source. (Provided for comparison use only).

PBL's 96-well ELISA kit will equip you with the most sensitive and consistent tool for accurately measuring Mouse IFN- β in serum and plasma matrices. Purified, recombinant mouse IFN- β expressed in mammalian cells is provided as the standard. The sensitivity provided by this kit will advance your mouse studies and help uncover insights into the pathology of diseases.

Figure 2. Precision and Recovery in CD-1 Mouse Serum

Concentration	2.5 pg/ml	10 pg/ml	50 pg/ml
Intra-Assay CV (n=24)	5.7%	5.5%	5.5%
Inter-Assay CV (n=9)	8.3%	6.3%	4.8%
Average Recovery (n=27)	98%	92%	94%

Figure 3. Comparison of ELISA Reactivity of Mammalian & *E. coli* Expressed Mu IFN- β



The versatility of mouse models lends their use in a number of disease and auto-immunity studies. Interferon β plays a pivotal role in the protective response to many infections and diseases due to its antiviral, anti-proliferative, anti-inflammatory and immune-modulatory activities, and is an important early product of TLR/RLR stimulation.

Product Information

Catalog Number	42410-1, -2
Description	VeriKine-HS™ Mouse Interferon Beta Serum ELISA
Size	1 x 96-well plate, 5 x 96-well plate
Compatibility	Serum, Plasma, Tissue Culture Media
Assay Range	0.94 - 60 pg/ml



For more information on this product, visit our website.