



# VeriKine™

## Mouse Interferon Alpha ELISA

Catalog No. 42120

- Provides global mouse IFN- $\alpha$  measurement in cell culture supernatant with the ability to detect all 14 mouse IFN- $\alpha$  subtypes
- Allows accurate mouse IFN- $\alpha$  measurement to 12.5 pg/ml
- Provides consistent performance and reproducible results with <10% inter- and intra-assay CV, as well as consistently low background
- Specific for mouse IFN- $\alpha$  with no cross-reactivity detected with rat IFN- $\alpha$ , mouse IFN- $\beta$  or  $\gamma$ , or human IFN- $\alpha$

**Figure 1.** VeriKine™ Mouse IFN- $\alpha$  ELISA subtype cross-reactivity

IFN- $\alpha$ Subtype	Results	Detection Limit (pg/ml)
42120 Standard	+++	12.5
IFN- $\alpha$ A	+++	12.5
IFN- $\alpha$ 14	+++	12.5
IFN- $\alpha$ 8	++	25
IFN- $\alpha$ 12	++	25
IFN- $\alpha$ 11	++	25
IFN- $\alpha$ B	++	50
IFN- $\alpha$ 1	+	50
IFN- $\alpha$ 5	+	50
IFN- $\alpha$ 6T	+	50
IFN- $\alpha$ 9	+	50
IFN- $\alpha$ 4	+	50
IFN- $\alpha$ 7	+	50
IFN- $\alpha$ 13	+	50
IFN- $\alpha$ 2	+	50

The VeriKine™ Mouse Interferon Alpha ELISA Kit has been developed to quantify levels of IFN- $\alpha$  in tissue culture media using a sandwich immunoassay. The kit contains all the reagents required to perform 1 x or 5 x 96-well plates.

PBL Assay Science (formerly PBL InterferonSource)  
131 Ethel Road West, Suite 6 | Piscataway, NJ 08854

**Phone:** (732) 777-9123

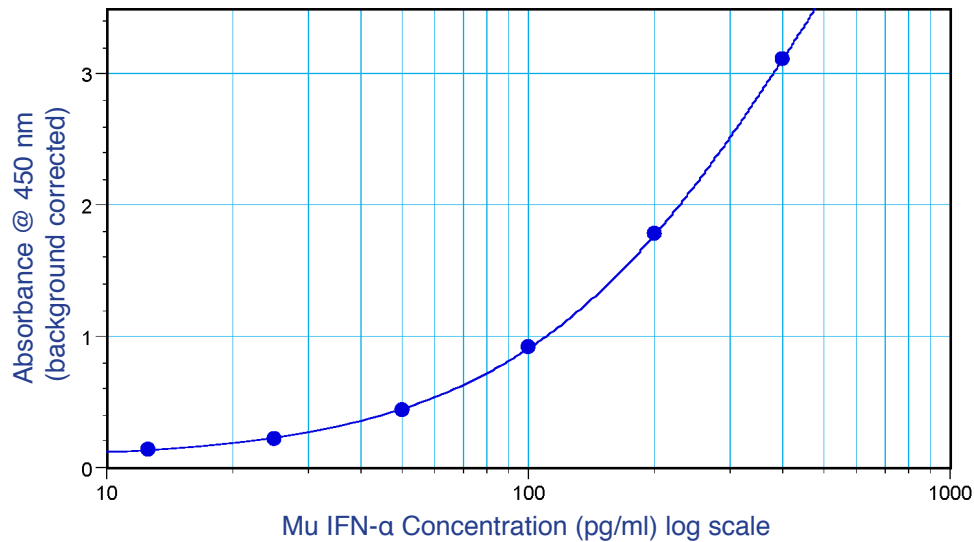
**Fax:** (732) 777-9141

**E-mail:** [info@pblassaysci.com](mailto:info@pblassaysci.com)

**Web:** [www.pblassaysci.com](http://www.pblassaysci.com)

In humans, Interferon Alpha (IFN- $\alpha$ ) protein therapeutics are used to treat viral diseases and cancer, and are often examined in mouse models of these diseases. Recent evidence has suggested that IFN- $\alpha$  proteins may play a role in the etiology of certain autoimmune diseases, and IFN- $\alpha$  therapeutic treatment may encompass a promising approach for other diseases.

**Figure 2. Representative Standard Curve of Cat. No. 42120**



### Selected Citations

1. E.W. Choi, *et al.* 2012. Reversal of serologic, immunologic, and histologic dysfunction in mice with systemic lupus erythematosus by long-term serial adipose tissue-derived mesenchymal stem cell transplantation. *Arthritis & Rheumatism* 64(1):243-253.
2. Lihoradovaa, O., *et al.* 2012. The Dominant-Negative Inhibition of Double-Stranded RNA-Dependent Protein Kinase PKR Increases the Efficacy of Rift Valley Fever Virus MP-12 Vaccine. *Journal of Virology* 86(14):7650-7661.

### Product Information

Catalog Number	42120-1, -2
Description	VeriKine™ Mouse Interferon Alpha ELISA
Size	1 x 96-well plate, 5 x 96-well plate
Compatibility	Cell Culture Supernatant
Assay Range	12.5 - 400 pg/ml

For more information and product availability, please contact our customer service department by telephone (732-777-9123) or e-mail ([info@pblsaysci.com](mailto:info@pblsaysci.com)).



For more information, including application notes, technical notes and product citations, visit our website.