



# VeriKine™

## Mouse Interferon Alpha ELISA

Catalog No. 42120

- Accurate mouse IFN- $\alpha$  measurement to 12.5 pg/ml in tissue culture media
- Ability to detect all 14 mouse IFN- $\alpha$  subtypes
- Consistent performance and reproducible results with < 10% inter- and intra-assay CVs
- Specific for mouse IFN- $\alpha$  with no cross-reactivity detected against 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

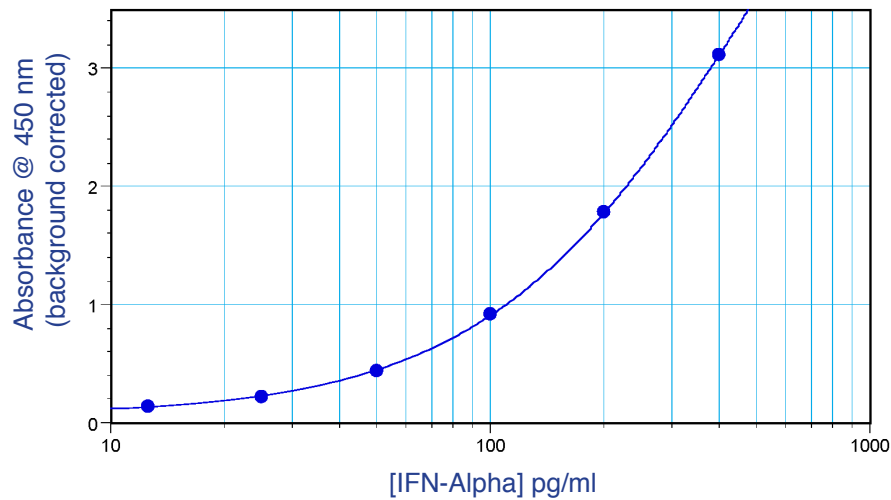
The VeriKine™ Mouse Interferon Alpha ELISA Kit has been developed to quantify levels of IFN- $\alpha$  in tissue culture media (TCM) by sandwich immunoassay. The kit contains all components required to perform 1 or 5 x 96-well plates.

**pbl** assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854

**T** (732) 777-9123 | **F** (732) 777-9141 | **E** info@pblassaysci.com | **W** www.pblassaysci.com

**Figure 2.** Representative Standard Curve of 42120 Mouse IFN Alpha



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.

### Product Information

Catalog Number	42120
Description	VeriKine™ Mouse Interferon Alpha ELISA
Size(s)	1 x 96-well plate (42120-1) 5 x 96-well plate (42120-2)
Compatibility	Tissue Culture Media (TCM)
Assay Range	12.5 - 400 pg/ml

**Visit our website for more information:**

<https://www.pblabsci.com/assay-kits/mouse-ifn-alpha-elisa-kit-tcm-42120>