

Certificate of Analysis

Human Interferon Alpha Sampler Set

Catalog No: 11002-1

Lot No:

Size: ~ 1 x 10⁵ units/vial

Introduction: While human interferon alpha A (Hu-IFN- A) has worked reasonably well in most applications, there are twelve identified individual species in the human interferon- family, each with vastly different properties. Some have proven to be ten and twenty times more effective than Hu-IFN- A in protecting a cell against specific viruses (see graph). Nonetheless, ninety-nine percent of the human interferon- sold for research is still IFN- A (IFN- 2a) or the closely related IFN- 2 (IFN- 2b).

The Human Interferon- Sampler can help you find the right alpha interferon for your research. The Sampler contains ~ 100,000 units of each of the species. By testing and selecting specific alpha interferons to meet your research needs, not only could you see improved effectiveness and reduce the amount of material required, but you could increase your understanding of how interferons function.

Product Information

Volume: ml

Activity: Refer to table for the activity (MDBK/VSV) of individual subtypes

Specific Activity: Refer to table for the specific activity (MDBK/VSV) of individual subtypes

Buffer: Phosphate buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

Endotoxin: < 1 EU/ g

Purity: ~ 95%

Purification Method: Purified from *E. coli* by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography

Source: cDNA obtained from human leukocyte mRNA expressed in *E. coli*

Assay Used to Measure Bioactivity: Interferon was titrated with the use of the cytopathic effect inhibition assay as listed

Bovine (MDBK/VSV) . performed as described [Rubinstein, *et al.* (1981) *J. Virol.* 37(2):755]. The EC₅₀ for interferon in this assay is ~5 U/ml. Lot Activity was derived from multiple determinations in the above assay.

Human (A549/EMCV) . performed as described [Budd, *et al.* (1985) *Canc. Chem. Pharm.* 12:39]. The EC₅₀ for interferon in this assay is ~1 U/ml.

The units are determined by use of a Human IFN Alpha A (Hu-IFN- A [2a]) laboratory standard calibrated to the international reference standard for Human Interferon Alpha A (Hu-IFN- A [2a]) provided by the National Institutes of

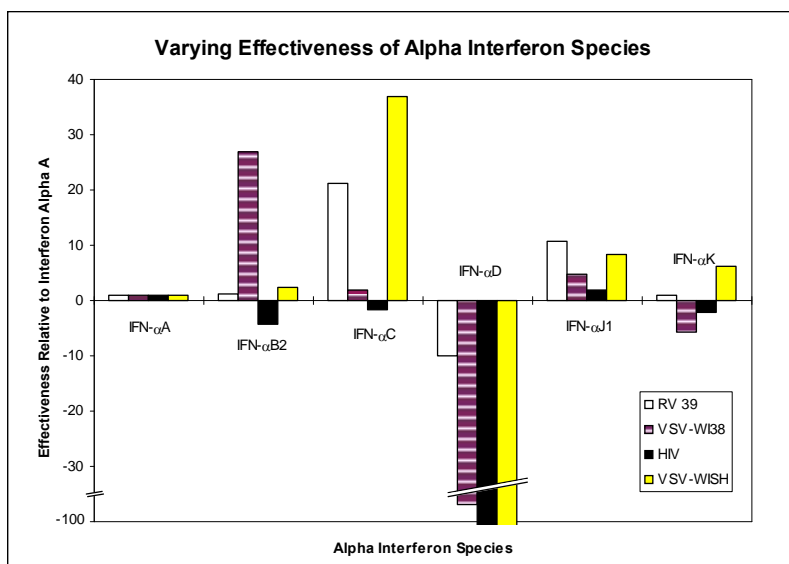


Figure 1. Varying Effectiveness of Alpha Interferon Species.
 1. Sperber, S., *et al.* (1992) *J. Interferon Research* 12:363-368.
 2. Sperber, S., *et al.* (1993) *Antiviral Research* 22:121-129.

Health [Meager, *et al.* (2001) *J. Immunol. Methods.* 257(1-2):17-33]. Please note that IFN assays vary between labs and assay systems [Meager, *et al.* (2001) *J. Immunol. Meth.* 257:17. Meager and Das (2005) *J. Immunol. Meth.* 306:1].

References: For more information on the various alleles of IFN alpha 2a, 2b, 4a, 4b, 1 and D.

1. Gewart, DR., *et al.* (1995) *J. Interferon Cytokine Res.* 15(5):403.
2. Hussain, M., *et al.* (1997) *J. Interferon Cytokine Res.* 17(9):559.
3. Hussain, M., *et al.* (2000) *J. Interferon Cytokine Res.* 20:763.

Materials Supplied

Part No.	Human Interferon Subtype Description	Component Lot No.	Accession Number	Gene	Molecular Weight (kDa)	Theoretical pI	Reference Catalog No.*
1	Alpha A (2a)		V00549	IFNA2	19.2	5.99	11100
2	Alpha B2 (8)		X03125	IFNA8	19.5	5.18	11115
3	Alpha C (10)		X02961	IFNA10	19.4	5.7	11120
4	Alpha D (1 [Val(114)])		V00538	IFNA1	19.4	5.18	11125
5	Alpha F (21)		V00540	IFNA21	19.3	5.99	11130
6	Alpha G (5)		X02956	IFNA5	19.5	5.46	11135
7	Alpha H2 (14)		X02959	IFNA14	19.7	6.38	11145
8	Alpha I (17)		V00532	IFNA17	19.3	5.46	11150
9	Alpha J1 (7)		X02960	IFNA7	19.6	5.87	11160
10	Alpha K (6)		X02958	IFNA6	19.7	6.43	11165
11	Alpha 4b (4)		X02955	IFNA4	19.4	5.76	11180
12	Alpha WA (16)		X02957	IFNA16	19.3	5.96	11190

Part No.	Human Interferon Subtype Description	Component Lot No.	Concentration (Activity) (MDBK/VSV) (U/ml)	Concentration (Activity) after 1x Freeze-Thaw cycle (MDBK/VSV) (U/ml)+	Activity on A549 (human lung epithelial cells)/EMCV (U/ml)	Specific Activity on MDBK/VSV (U/mg)
1	Alpha A (2a)		x 10	x 10	x 10	x 10
2	Alpha B2 (8)		x 10	x 10	x 10	x 10
3	Alpha C (10)		x 10	x 10	x 10	x 10
4	Alpha D (1 [Val(114)])		x 10	x 10	x 10	x 10
5	Alpha F (21)		x 10	x 10	x 10	x 10
6	Alpha G (5)		x 10	x 10	x 10	x 10
7	Alpha H2 (14)		x 10	x 10	x 10	x 10
8	Alpha I (17)		x 10	x 10	x 10	x 10
9	Alpha J1 (7)		x 10	x 10	x 10	x 10
10	Alpha K (6)		x 10	x 10	x 10	x 10
11	Alpha 4b (4)		x 10	x 10	x 10	x 10
12	Alpha WA (16)		x 10	x 10	x 10	x 10

* The Reference Catalog Numbers are provided for ease of ordering the individual human alpha subtypes. Please use this information to order individual alpha subtypes. Additional background information on the different interferon alpha subtypes is available upon request.

+ See Storage Conditions/Comments.



Shipping Information

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, the product may be stored at -20°C for short-term use (m6 months). For long-term storage, we recommend storing the product at -70°C or below for retention of full activity. When thawing, the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA) or tissue culture media with serum. Dilution of material below 2×10^5 units/ml for freezing is not recommended. There may be losses in activities after one freeze-thaw cycle. One freeze-thaw cycle is equivalent to thawing an aliquot prepared from the material received. Refer to the table for activities on MDBK/VSV cells post one freeze-thaw cycle. Please note that IFN activities vary across labs and assay systems. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____

Date:

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.

Rev. 04