

Human Interferon Alpha Sampler Set

This unique product contains a set of 12 different recombinant human IFN- α species. Some of them have proven to be ten and twenty times more effective than human IFN- α A in protecting a cell against specific viruses (see table). It is a ideal tool for low price first screening experiments and allows quick decisions on further research directions.

Ordering Information

[Order Online »](#)

PBL-11002-1	1Set
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon- α , IFN- α

Contents

- IFN- α A (Prod. No. [PBL-11100](#))
- IFN- α B2 (Prod. No. [PBL-11115](#))
- IFN- α C (Prod. No. [PBL-11120](#))
- IFN- α D (Prod. No. [PBL-11125](#))
- IFN- α F (Prod. No. [PBL-11130](#))
- IFN- α G (Prod. No. [PBL-11135](#))
- IFN- α H2 (Prod. No. [PBL-11145](#))
- IFN- α I Prod. No. [PBL-11150](#))
- IFN- α J1 (Prod. No. [PBL-11160](#))
- IFN- α K (Prod. No. [PBL-11165](#))
- IFN- α 4b (Prod. No. [PBL-11180](#))
- IFN- α WA (Prod. No. [PBL-11190](#))

Endotoxin Content < 1 EU/ μ g

Purity \geq 95%

Purity Detail Purified from E. coli by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography

Quantity 1×10^5 U/vial.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com