

Mouse Interferon Alpha A

Citations: 5

[View Online »](#)

Ordering Information

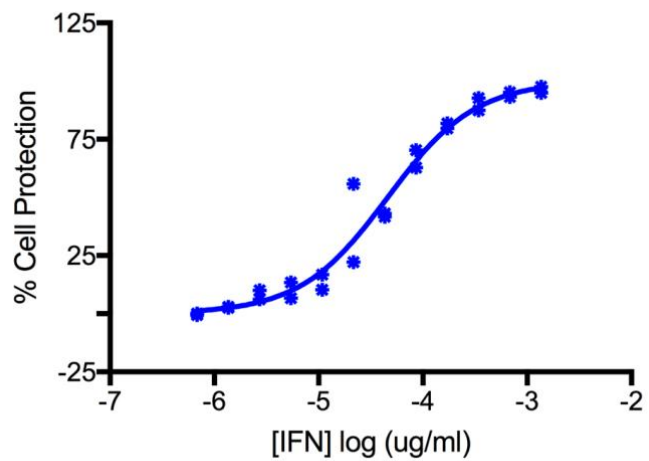
[Order Online »](#)

PBL-12100-1	0.1MU
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Mouse IFN Alpha A recombinant protein with activity > 1 x 10⁵ U/ml
- Distinct IFN-alpha subtype used extensively for research
- Greater than 95% purity



Antiviral Activities of Mouse IFN-Alpha A (12100) and Mouse IFN-Alpha 4 (12115) on Mouse L929 Cells infected with EMCV

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 A, IFN- α A
Endotoxin Content	< 1 EU/ μ g
Formulation	Liquid. In PBS containing 0.1% BSA.
MW	~19.3kDa
Purity	\geq 95%
Purity Detail	A combination of ion exchange, hydrophobic interaction and size exclusion chromatography.
Quantity	0.1ml
Source	Produced in <i>E. coli</i> .



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