

Human Interferon Alpha A (Alpha 2a)

Citations: 4

[View Online »](#)

Ordering Information

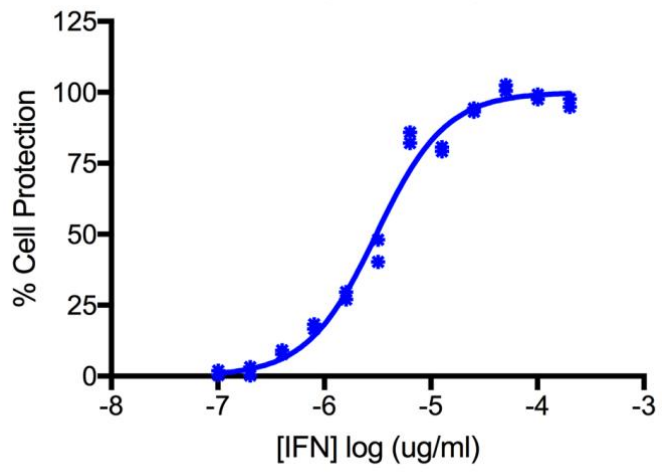
[Order Online »](#)

PBL-11100-1	5x1MU
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

- Human IFN Alpha 2a recombinant protein with high activity
- Less than 1 EU/μg endotoxin
- 95% or greater purity



Titration of Human IFN-Alpha 2a in A549 Cells

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 Aa, IFN- α 2a, Interferon- α 2a

Endotoxin Content < 1 EU/ μ g

Formulation Liquid. In PBS containing 0.1% BSA.

MW ~19.2kDa

Purity \geq 95%

Purity Detail A combination of ion exchange, hydrophobic interaction and size exclusion chromatography.

Quantity 0.1ml

Source Produced in *E. coli*.



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