

Universal Type I Interferon

IFN- α hybrid constructed from recombinant human IFN- α A and IFN- α D. Active on almost every mammalian cell, substituting type I IFN- α , IFN- β , IFN- ω and IFN- τ .

Citations: 5

[View Online »](#)

Ordering Information

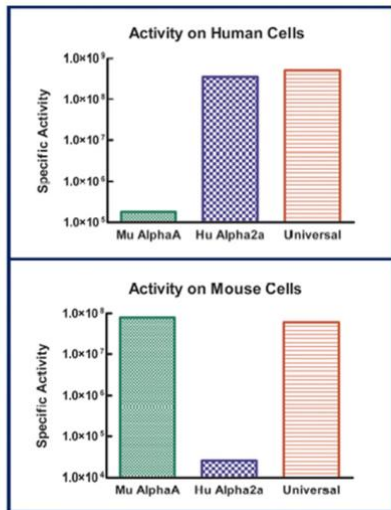
[Order Online »](#)

PBL-11200-1	0.1MU
PBL-11200-2	1MU

Manuals, SDS & CofA

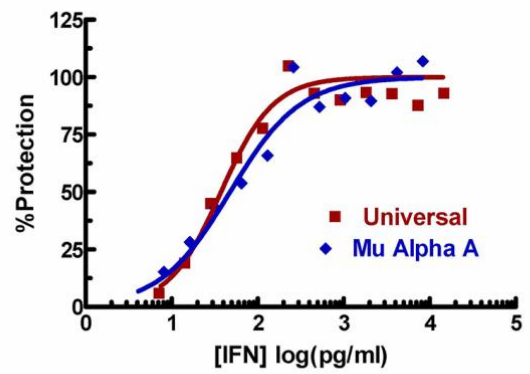
[View Online »](#)

- Recombinant Human IFN-Alpha protein hybrid exhibiting activity across multiple species
- Crucial for cross-species comparative studies in virology, immune modulation and more
- Essential to *in vivo* or *in vitro* studies where the species-relevant IFN alpha is unavailable or performs inconsistently



Activity Comparison of Universal™ Type I IFN, Human IFN-Alpha 2a and Mouse IFN-Alpha A on Human and Mouse Cell Lines

Comparison of Universal IFN and Mouse Alpha A Antiviral Activity on Mouse Cells



Antiviral Activities of Universal™ Type I IFN (11200) and Mouse IFN-Alpha A 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 Universal Type I Interferon, IFN- α A/D (BgIII), Interferon- α A/D (BgIII)

Endotoxin Content < 1 EU/ μ g

Formulation Liquid. In PBS containing 0.1% BSA.

MW ~19.4kDa

Purity \geq 95%

Purity Detail Purified from *E. coli* by 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