

C12R Mouse IFN Alpha Decoy Receptor, neutralizing

C12R is an interferon binding/blocking protein produced by many strains of Mousepox (Ectromelia) virus. This protein blocks activity of mouse IFN Alpha but not Beta. It is an essential protein for viral infection.

Ordering Information

[Order Online »](#)

PBL-12185-1	10µg
PBL-12185-2	100µg

Manuals, SDS & CofA

[View Online »](#)

- Viral C12R IFN Antagonist selectively neutralizes **all** mouse IFN alpha subtypes
- No cross-reactivity with mouse IFN beta
- Also neutralizes all human IFN alpha subtypes, human IFN beta, human IFN omega, rat IFN alpha 1, and mouse limitin
- Suitable for Western Blot

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Biological Activity Measured by neutralization of interferon in cytopathic effect inhibition assay.

Endotoxin Content < 1 EU/μg

Formulation Phosphate-buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

Purity > 95%

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

Specificity Neutralizes all mouse IFN alpha subtypes, all human IFN alpha subtypes, human IFN beta, human IFN omega, rat IFN alpha 1, and mouse limitin.
Does not neutralize mouse IFN beta, human IFN gamma, human IFN lambda 1, human IFN lambda 2, human IFN lambda 3, or rat IFN alpha 14 when tested at ≥ 100X the typical neutralizing dose.



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