

Porcine Interferon Alpha

Ordering Information

[Order Online »](#)

PBL-17105-1	0.1MU
-------------	-------

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

Endotoxin Content < 1 EU/ μ g

Formulation Liquid. In PBS containing 0.1% BSA.

Purity \geq 95%

Purity Detail Purified from mammalian tissue culture media by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography.

Quantity 0.1ml

Source Produced in mammalian cells.



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