

Mouse Interferon Beta, mammalian, carrier-free

- Authentic sequence of mouse interferon beta protein with no tag

- 10 fold greater specific activity than standard protein

Ordering Information

[Order Online »](#)

PBL-12410-1	1MU
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	After receipt, this product should be kept at -70°C or below for retention of full activity. Thaw product vial by incubation in cold tap water until just thawed – the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA).
Handling	After opening, prepare aliquots and store at -80°C. Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Interferon-β, IFN-β
Appearance	Liquid
Endotoxin Content	< 1 EU/μg
Formulation	Liquid. 20 mM Sodium Phosphate, pH 6.4; 300 mM NaCl; 5% Glycerol.
MW	~19.6kDa
Purity	≥95%
Purity Detail	Purified from mammalian tissue culture media by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography.
Quantity	≥ 1 x 10 ⁶ units
Source	Gene was obtained from mouse DNA expressed in mammalian cells



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)