

Mouse Interferon Beta, mammalian

- Authentic sequence of mouse interferon beta protein
- 10 fold greater specific activity than E. coli derived protein

Ordering Information

[Order Online »](#)

PBL-12405-1	0.1MU
-------------	-------

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
Biological Activity	<p>Assay Used to Measure Bioactivity: Interferon was titrated with the use of the cytopathic effect inhibition assay as listed. Mouse (L929/EMCV)-performed as described [Rubinstein et al. (1981) J. Virol. 37 (2):755]. The EC50 for interferon in this assay is ~ 2.5 U/ml. Lot Activity was derived from multiple determinations in the above assay.</p> <p>The units are determined by use of a Mu IFN-β laboratory standard calibrated to the international reference standard for mouse Interferon Beta provided by the National Institutes of Health [Meager et al. (2001) J. Immunol Methods. 257(1-2):17-33]. Please note that IFN assays vary between labs and assay systems [Meager et al (2001). J. Immunol. Meth. 257:17. Meager and Das (2005) J. Immunol. Meth. 306:1].</p>
Endotoxin Content	< 1 EU/ μ g
Formulation	Liquid. 20 mM Sodium Phosphate, pH 6.4; 300 mM NaCl; 5% Glycerol, 0.1% bovine serum albumin (BSA).
MW	~19.6kDa
Purity	\geq 95%
Purity Detail	Purified from mammalian tissue culture media by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography.

Quantity

$\geq 1 \times 10^5$ 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)