

# Mouse Interferon Beta, carrier-free

## Ordering Information

[Order Online »](#)

|             |       |
|-------------|-------|
| PBL-12401-1 | 0.1MU |
|-------------|-------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                          |  |
|--------------------------|--|
| <b>Use/Stability</b>     | Thaw product vial by incubation in cold tap water until just thawed. 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). |
| <b>Handling</b>          | After opening, prepare aliquots and store at -80°C. Avoid freeze/thaw cycles.  |
| <b>Long Term Storage</b> | -80°C  |
| <b>Shipping</b>          | Dry Ice  |

**Regulatory Status** RUO - Research Use Only

## Product Details

|                          |   |
|--------------------------|---|
| <b>Alternative Name</b>  | Interferon- $\beta$ , IFN- $\beta$  |
| <b>Endotoxin Content</b> | < 1 EU/ $\mu$ g   |
| <b>Formulation</b>       | Liquid. In PBS.   |
| <b>MW</b>                | ~19.6kDa  |
| <b>Purity</b>            | $\geq$ 95%  |
| <b>Purity Detail</b>     | A combination of ion exchange, hydrophobic interaction and size exclusion chromatography. |
| <b>Source</b>            | Produced in <i>E. coli</i> .  |



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)