

# Anti-Human Interferon Beta, Clone MMHB-12 (MAb)

Neutralizes human interferon beta; binds to human interferon beta with high affinity

## Ordering Information

[Order Online »](#)

|             |      |
|-------------|------|
| PBL-21450-1 | 50µg |
|-------------|------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |   |
|-------------------|---|
| Handling          | After opening, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles. |
| Long Term Storage | -20°C   |
| Shipping          | Dry Ice   |

**Regulatory Status** RUO - Research Use Only

## Product Details

|                    |  |
|--------------------|--|
| Alternative Name   | IFN- $\beta$ , Interferon- $\beta$         |
| Clone              | MMHB-12                                    |
| Formulation        | Liquid. In PBS containing 0.1% BSA.        |
| Host               | Mouse                                      |
| Immunogen          | Human IFN- $\beta$ (interferon- $\beta$ ). |
| Isotype            | IgG1                                       |
| Species Reactivity | Human                                      |



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)