

# Anti-Human Interferon Gamma Receptor Chain 2, Clone MMHGR-2 (MAb)

Binds to but does not neutralize human interferon gamma receptor

## Ordering Information

[Order Online »](#)

|             |       |
|-------------|-------|
| PBL-21585-1 | 250µg |
|-------------|-------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

|                           |   |
|---------------------------|---|
| <b>Alternative Name</b>   | IFN- $\gamma$ Receptor Chain 2, Interferon- $\gamma$ Receptor Chain 2 |
| <b>Clone</b>              | MMHGR-2   |
| <b>Formulation</b>        | Liquid. In PBS containing 0.1% BSA.                                   |
| <b>Host</b>               | Mouse   |
| <b>Immunogen</b>          | Human IFN- $\gamma$ (interferon- $\gamma$ ) receptor chain 2.         |
| <b>Isotype</b>            | IgG1  |
| <b>Species Reactivity</b> | 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)