

Anti-Human Interferon Alpha, Clone MMHA- 11 (MAb)

Neutralizes human interferon alpha; binds to human interferon alpha

Ordering Information

[Order Online »](#)

| | |
|-------------|------|
| PBL-21112-1 | 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 | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------|--------------------------------------|
| Alternative Name | IFN- α , Interferon- α |
| Application | ELISA |
| Clone | MMHA-11 |
| Formulation | Liquid. In PBS containing 0.1% BSA. |
| Host | Mouse |
| Immunogen | Human leukocyte interferon. |
| Isotype | IgG1 |
| Species Reactivity | Human |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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