

# Anti-Mouse Interferon Alpha, Clone RMMA-1 (MAb)

Neutralizes mouse interferon alpha; binds to mouse interferon alpha

## Ordering Information

[Order Online »](#)

PBL-22100-3	FITC - 25µg
PBL-22100-1	250µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	PBL-22100-1 (Unlabelled) should be shipped on dry ice. Storage -20°C. PBL-22100-3 (FITC-labelled) should be shipped on blue ice. Storage +4°C. Protect from prolonged exposure to light.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	IFN- $\alpha$ , Interferon- $\alpha$
<b>Application</b>	ELISA
<b>Clone</b>	RMMA-1
<b>Formulation</b>	Liquid.
<b>Host</b>	Rat
<b>Immunogen</b>	Mouse IFN- $\alpha$ (interferon- $\alpha$ ).
<b>Isotype</b>	IgG1
<b>Species Reactivity</b>	Mouse



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)