

# Mouse Monoclonal Antibody against Human Interferon Alpha/Beta Receptor

## 1, Clone MMHAR-3

Detects human interferon alpha/beta R1

**(MAb)**

Ordering Information

[Order Online »](#)

PBL-21370-1	50µg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Handling</b>	For long term storage keep unconjugated antibody at -20°C. Avoid freeze/thaw cycles. Keep conjugated formats at +4°C. Do not freeze. Protect from light.
<b>Short Term Storage</b>	-20°C
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	IFN- $\alpha/\beta$ Receptor Chain 1, Interferon- $\alpha/\beta$ Receptor Chain 1
<b>Application</b>	Flow Cytometry
<b>Clone</b>	MMHAR-3
<b>Formulation</b>	Phosphate – buffered saline (PBS) containing 1 .5% BSA and 0.01 % Kathon as a preservative, pH 7.4
<b>Host</b>	Mouse
<b>Immunogen</b>	Human Interferon alpha/beta receptor 1
<b>Isotype</b>	IgG1



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-fr@enzolifesciences.com](mailto:info-fr@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)