

# IL-6 receptor Recombinant monoclonal antibody (rhPM-1) (Human IgG1κ)

The IL-6 receptor is a protein complex consisting of an IL-6 receptor subunit (IL6R) and interleukin 6 signal transducer Glycoprotein 130. Interleukin 6 (IL6) is a potent pleiotropic cytokine that regulates cell growth and differentiation and plays an important role in immune response.

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

ENZ-ABS551-0200	200µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Interleukin-6 receptor, IL6R, MRA
Application	ELISA
Clone	rhPM-1 (Tocilizumab)
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Human
Immunogen	This antibody was prepared by CDR grafting of mouse PM-1 antibody onto a human REI and NEW framework regions to create rhPM-1.
Isotype	IgG1k
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	Recognizes both the membrane bound and soluble forms of human IL-6R.
UniProt ID	P08887

### Worry-free Guarantee



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)

### This antibody is covered by our **Worry-Free Guarantee**.