IL-6 receptor Recombinant monoclonal antibody (rhPM-1) (Rabbit IgGκ)

The IL6 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-ABS550-0200

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 Rabbit

Immunogen The original antibody was generated by CDR grafting of

mouse PM-1 antibody onto human REI and NEW

framework regions to create rhPM-1.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

Specificity Recognizes both the membrane bound and soluble forms

of human IL-6R.

Technical Info / Product NotesThis chimeric rabbit antibody was made using the variable

domain sequences of the original Human IgG1 format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID A0N0L5

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: August 5, 2024



info-

eu@enzolifesciences.com