TNF-R2 monoclonal antibody (R00N6)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

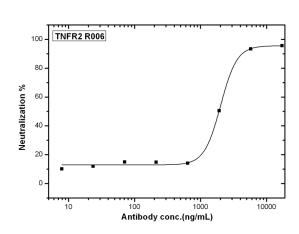
Order Online »

ENZ-ABS950-0200

200µg

Manuals, SDS & CofA

View Online »



TNFR2/TNFRSF1B-mediated inhibition of cytotoxicity was Neutralized by Human TNFR2 Antibody. Recombinant Human TNFR2/TNFRSF1B inhibits Recombinant Human TNFa induced cytotoxicity in the L-929 mouse fibroblast cell line. Inhibition of Recombinant Human TNFa (0.25 ng/mL) activity elicited by Recombinant Human TNFR2/TNFRSF1B (1 μ g/mL) is neutralized by increasing concentrations of Human TNFR2/TNFRSF1B Monoclonal Antibody. The IC50 is typically 1-3 μ g/mL in the presence of the metabolic inhibitor actinomycin D (1 μ g/mL).

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD120b, p75, TNF-R 80kDa, TNFRSF 1B

Application Neutralization

Clone R00N6

Formulation Liquid. 0.2 µm filtered solution in Histidine and Arginine

buffer containing 120mM NaCl, 0.02% Tween 80, pH 6.0.

Preservative free.

Host Rabbit

Immunogen Recombinant human TNFR2 protein.

lgG lgG

Purity Detail Protein A purified.

Species Reactivity Human

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

_

Last modified: September 17, 2024

