TNF-R1 monoclonal antibody (R111)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

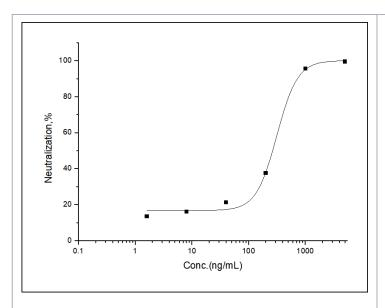
Order Online »

ENZ-ABS964-0200

200µg

Manuals, SDS & CofA

View Online »



Recombinant Human TNFR1/TNFRSF1A inhibits Recombinant Human TNFa induced cytotoxicity in the L-929 mouse fibroblast cell line. Inhibition of Recombinant Human TNFa (0.2 ng/mL) activity elicited by Recombinant Human TNFR1/TNFRSF1A (0.3 μ g/mL) is neutralized by increasing concentrations of Human TNFR1/TNFRSF1A Monoclonal Antibody. The IC50 is typically 0.2-1 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 p55, CD120a, TNF-R 60kDa, TNFRSF 1A

Application Neutralization

Clone R111

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 TNFR1 / CD120a / TNFRSF1A

protein.

Isotype IgG

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



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com