

TNF-R2 monoclonal antibody (R00N6)

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

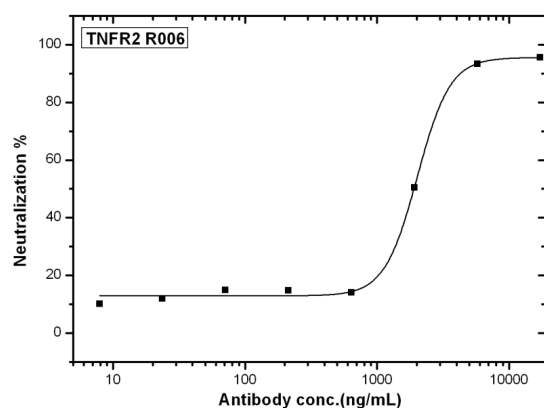
Ordering Information

[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 TNF α induced cytotoxicity in the L-929 mouse fibroblast cell line. Inhibition of Recombinant Human TNF α (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.

Isotype IgG

Purity Detail Protein A purified.

Species Reactivity Human

Worry-free Guarantee This antibody is covered by our [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)