

TRAIL-R2 polyclonal antibody

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

Citations: 20

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TRAIL receptor 2, DR5, Death receptor 5, KILLER, CD262, Trick2, Ly98, TNFRSF10B, TNF-related apoptosis-inducing ligand receptor 2, Tumor necrosis factor receptor superfamily member 10B
Application	Flow Cytometry, ICC, IHC (PS), WB
Application Notes	Detects a band of ~52kDa by Western blot. Cited applications include Immunoprecipitation.
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Gene/Protein Identifier	8795 (Entrez GeneID)
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 68-94 (A68PQQKRSSPSEGLCPPGHHISEDGRDC94) of the extracellular domain of human TRAIL-R2.
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Western Blot or Immunocytochemistry on cell lines HEK 293 or Jurkat.

Recommendation Dilutions/Conditions

Immunocytochemistry (1:300)Immunohistochemistry (paraffin sections)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Human, Mouse

UniProt ID

O14763

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com