

TRAIL-R1 polyclonal antibody

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

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TRAIL receptor 1, Apo-2, CD261, DR4, Death receptor 4, TNFRSF 10A, TNF-related apoptosis-inducing ligand receptor 1, Tumor necrosis factor receptor superfamily member 10A
Application	WB
Application Notes	Detects a band of ~57kDa by Western blot.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
GenBank ID	U90875
Gene/Protein Identifier	8797 (Entrez GeneID)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 427-445 of human TRAIL-R1.
Species Reactivity	Human, Mouse, Rat
UniProt ID	O00220
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](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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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