

# TRAIL-R1 polyclonal antibody

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

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-SA225-0100	100µg
----------------	-------

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 <a href="#">Worry-Free Guarantee</a>

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-be@enzolifesciences.com](mailto:info-be@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-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)