

# TRAIL (human) monoclonal antibody (RIK-2)

Induces apoptosis in tumorigenic and transformed cell lines, involved in T cell mediated cytotoxicity, TRAIL ligand of cognate receptors induces NF-kappa-B signaling in a variety of cell types. TRAIL is a 30kDa protein known as TNF-related apoptosis inducing ligand, CD253, TNFSF10, or Apo-2L. It can be expressed as a cell surface (30kDa) and soluble ligand (19-20kDa) produced by megakaryocytes and platelets, monocytes, neutrophils, NK cells and activated T cells. This antigen is also expressed on tumor cells and transformed cell lines. TRAIL has been reported to induce apoptosis in tumor and transformed cell lines by a caspase-dependent process. This antibody is useful for flow cytometric staining and blocking studies.

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

## Ordering Information

[Order Online »](#)

ALX-804-907-0100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Alternative Name	Apo-2L, TNFSF 10, CD253
Application	Flow Cytometry
Clone	RIK-2
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.
GenBank ID	U57059
Gene/Protein Identifier	8743 (Entrez GeneID)
Host	Mouse
Immunogen	Human TRAIL transfected mouse cell line.
Isotype	IgG1
Purity Detail	Affinity purified.
Species Reactivity	Human
UniProt ID	P50591
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-  
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)