

TRAIL-R3 (human) polyclonal antibody

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

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-744-C200	200µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

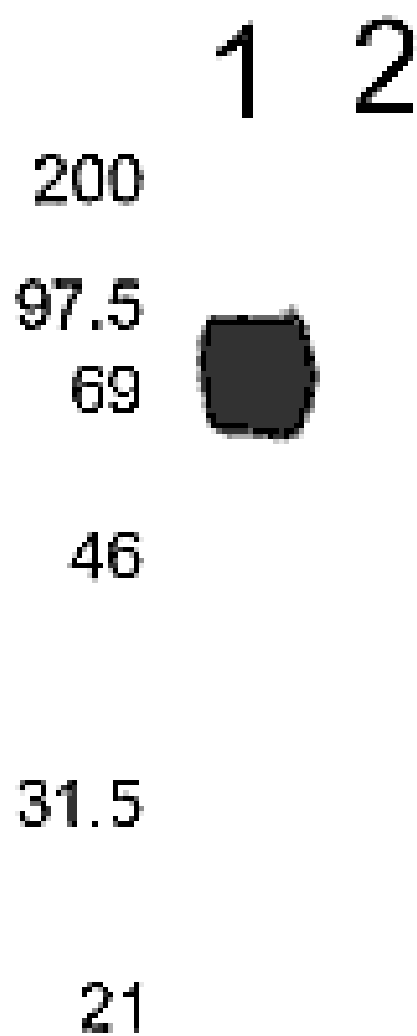


Figure: Specificity of PAb to TRAIL-R3. Lane 1: TRAIL-R3:Fc. Lane 2: TRAIL-R2. ALX-210-744 recognizes an approx. 66kDa band of the expected size of the recombinant TRAIL-R3:Fc protein in lane 1 and shows no detectable cross-reactivity to TRAIL-R2:Fc in lane 2.

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.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TRAIL receptor 3, DcR1, Death Receptor 1, TRID, CD263, TNFRSF 10C, TNF-related apoptosis-inducing ligand receptor 3, Tumor necrosis factor receptor superfamily member 10C
Application	Flow Cytometry, ICC, IHC (PS), WB
Application Notes	Detects a band of ~33kDa by Western blot.
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Gene/Protein Identifier	8794 (Entrez GeneID)
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 33-63 (E33VPQQTVAPQQQRHSFKGEECPAGSHRSEHT63) of the extracellular domain of human TRAIL-R3.
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Western blot or Immunocytochemistry on cell lines HEK 293 or Jurkat.
Recommendation	Flow Cytometry (1µl/106 cells/100µl)Immunocytochemistry (1:300),Western Blot
Dilutions/Conditions	(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



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

014798

European Sales Office
ENZO LIFE SCIENCES
(EU) 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

Worry-free Guarantee

This antibody is covered by our **Worry-Free Guarantee**.

For local distributors and detailed product information visit us online: www.enzo.com All product names, logos, and brands are the property of their respective owners.