

# TRAIL (human) polyclonal antibody

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

Citations: 5

[View Online »](#)

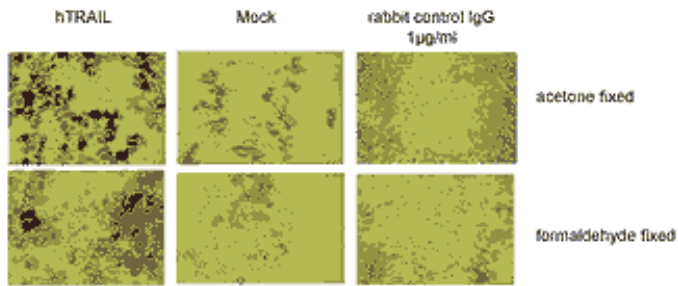
## Ordering Information

[Order Online »](#)

ALX-210-732-R100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



**Figure 3:** Immunostaining of HEK 293T cells transfected with a expression plasmid coding for human TRAIL (left panels), or mock transfected (right panel).

**Method:** 3 days after transfection of cells as indicated, cells were fixed with 4% formaldehyde or acetone for 5 min. at RT. Slides were blocked with IgG, and incubated for 1 hour with PAb to TRAIL (Prod. No. [ALX-210-732](#)) at a dilution of 1:100 or control rabbit IgG in PBS containing 1% BSA. After washes with PBS, samples were incubated with a biotinylated secondary antibody to rabbit IgG for 1 hour, washed in PBS and revealed with StreptABComplex/HRP (Vector) and AEC. """, TRAIL (human) polyclonal antibody Immunocytochemistry, TRAIL (human) polyclonal antibody Immunocytochemistry, yes ALX-210-732, 210-

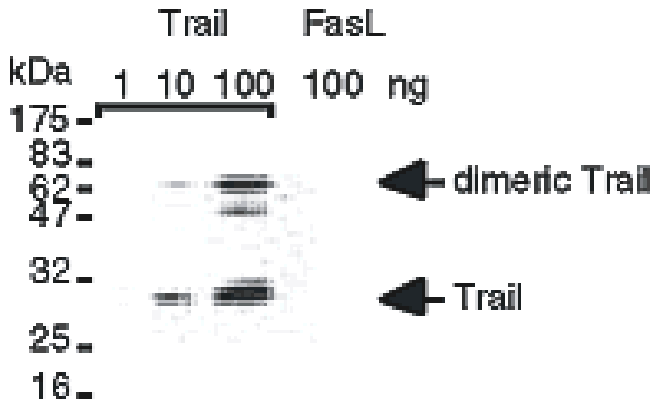
732\_immunogen.gif, [http://static.enzolifesciences.com/fileadmin/files/image/210-](http://static.enzolifesciences.com/fileadmin/files/image/210-732_immunogen.gif)

732\_immunogen.gif, **Figure 1:** Schematic structure of the immunogen., TRAIL (human) polyclonal antibody Schematic structure of immunogen, TRAIL (human) polyclonal antibody Schematic structure of immunogen, yes

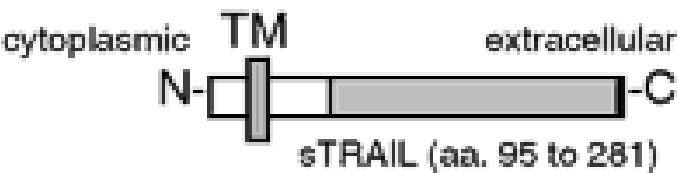
ALX-210-732, 210-

732.gif, [http://static.enzolifesciences.com/fileadmin/files/image/210-](http://static.enzolifesciences.com/fileadmin/files/image/210-732.gif)

732.gif, **Figure 2:** Detection of recombinant human TRAIL (Prod. No. [ALX-522-003](#)) but not of recombinant human FasL (Prod. No. ALX-522-001). **Method:** Recombinant soluble TRAIL



**Figure 2:** Detection of recombinant human TRAIL (Prod. No. ALX-522-003) but not of recombinant human FasL (Prod. No. ALX-522-001). **Method:** Recombinant soluble TRAIL, and recombinant soluble FasL as negative control, were resolved by SDS-PAGE under reducing conditions, transferred to nitrocellulose and detected with PAb to TRAIL (human) (Prod. No. ALX-210-732) at a dilution of 1:500. Proteins were visualized using a peroxidase-conjugated antibody to rabbit IgG and a chemiluminescence detection system.



**Figure 1:** Schematic structure of the immunogen.

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
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, ICC, IHC (FS), WB
Crossreactivity	Does not cross-react with mouse TRAIL.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
GenBank ID	U57059
Gene/Protein Identifier	8743 (Entrez GeneID)
Host	Rabbit
Immunogen	Recombinant human soluble TRAIL (aa 95-281).
Purity Detail	Protein A-affinity purified.
Source	Purified from serum.
Species Reactivity	Human
Specificity	Recognizes soluble TRAIL.
UniProt ID	P50591



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 940.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(EU) LTD  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
(EU) LTD  
Phone: +32 23 966 0426  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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

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

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

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