

TRAIL (human) polyclonal antibody

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

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

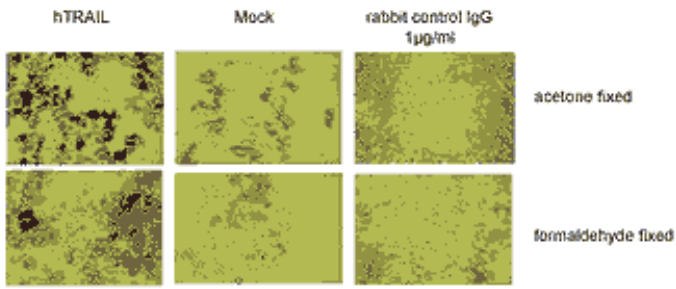


Figure 3: Immunostaining of HEK 293T cells transfected with a expression plasmid coding for human TRAIL (left panels), or mock transfected (right panel).

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.

Immunocytochemistry, TRAIL (human) polyclonal antibody Immunocytochemistry, yes ALX-210-732,210-732_immunogen.gif,http://static.enzolifesciences.com/fileadmin/files/image/210-732_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/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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.

ALX-210-732,210-732.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-732.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, 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 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.

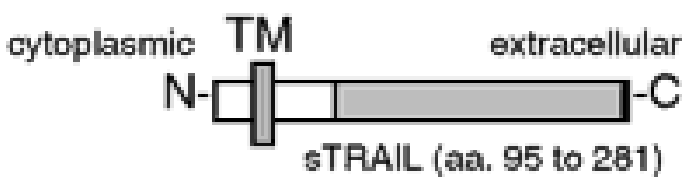


Figure 1: Schematic structure of the immunogen.

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
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-
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)