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 »

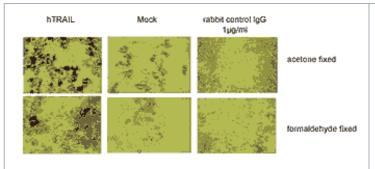


Figure 3: Immunostaining of HEK 293T cells transfected with a expression plasmid coding for human TRAIL (left panels), or mock transfected (right panel).

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-

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

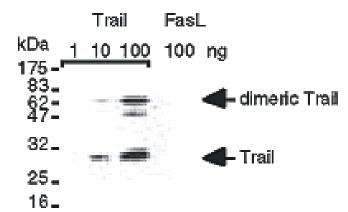


Figure 2: Detection of recombinant human TRAIL (Prod. No. ALX-522-003) but not of recombinant human FasL (Prod. No. ALX-522-001). Method: Recombinant soluble TRAIL, and recombinant soluble FasL as negative control, were resolved by SDS-PAGE under reducing conditions, transferred to nitrocellulose and 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.

cytoplasmic TM extracellular
NsTRAIL (aa. 95 to 281)

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 GenelD)

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



info-

eu@enzolifesciences.com