# TRAIL (soluble) (human), (recombinant) (Histag)

TRAIL, also called Apo2L, is a member of the tumor necrosis factor family (TNF). In contrast to other family members such as TNFa/TNFR1 and FasL/Fas, TRAIL and its receptors DR4 and DR5 are widely expressed in many human tissues. TRAIL activates rapid apoptosis in many transformed cell lines but not in normal tissues, even though TRAIL and its receptors DR4 and DR5 are expressed on both cell types. A glycophospholipid-anchored cell surface protein designated decoy receptor 1 (DcR1 or TRID) acts as a decoy to inhibit TRAIL signaling. Thus, TRAIL and its signaling pathways may have direct applications in selective tumor cell death and cancer research.

Citations: 48

View Online »

# **Ordering Information**

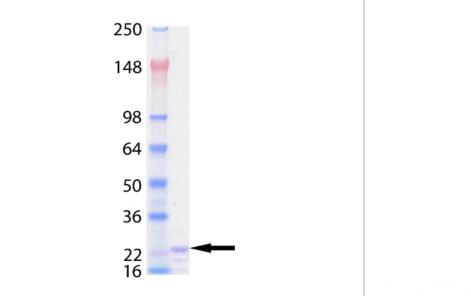
**Order Online** »

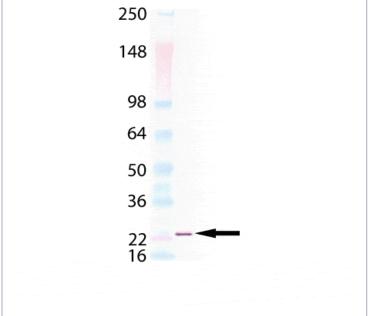
BML-SE721-0100

100µg

Manuals, SDS & CofA

**View Online** »





SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 1µg TRAIL (soluble) (human), (recombinant) (His-tag) (Prod. No. BML-SE721).

Western Blot Analysis: Lane 1: MW marker, Lane 2: 50 ng of TRAIL (soluble) (human), (recombinant) (His-tag) probed with Anti-6X His tag® antibody.

## **Handling & Storage**

**Use/Stability** When stored at -80°C, this protein is stable at the concentration supplied, in its current

storage buffer. Procedures such as dilution of the protein, followed by refreezing, could

lead to loss of activity.

**Handling** After initial defrost, aliquot product into individual tubes and refreeze at -70°C. Avoid

freeze/thaw cycles.

**Long Term Storage** -80°C

Shipping Dry Ice

### Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Apo-2L, TNFSF 10, CD253

Biological Activity Induces apoptosis in U937 cells at 1µg/ml, within 5 hours.

At 0.5-2.0µg/ml, induces apoptosis in various thyroid carcinoma lines within 18 hours. User must determine optimal concentrations and time course for induction. Please note that this protein does not require a cross-

linking enhancer to exhibit this potent activity.

**Endotoxin Content** <0.1 EU/μg TRAIL, determined using chromogenic

Limulus amebocyte lysate (LAL) test.

**Formulation** Liquid. In PBS, containing 10% glycerol and 0.1mM DTT.

**MW** ~23kDa

Purity ≥95% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Source Produced in E. coli. Human TRAIL (aa 114-281) is fused to

a His-tag.

Specificity Binds to human and mouse TRAIL receptors and human

osteoprotegerin (OPG).

UniProt ID P50591

Last modified: May 29, 2024



info-

eu@enzolifesciences.com