

# TRAIL (soluble) (mouse), (recombinant)

Ligand for TRAIL receptor (DR4, DR5).

Citations: 12

[View Online »](#)

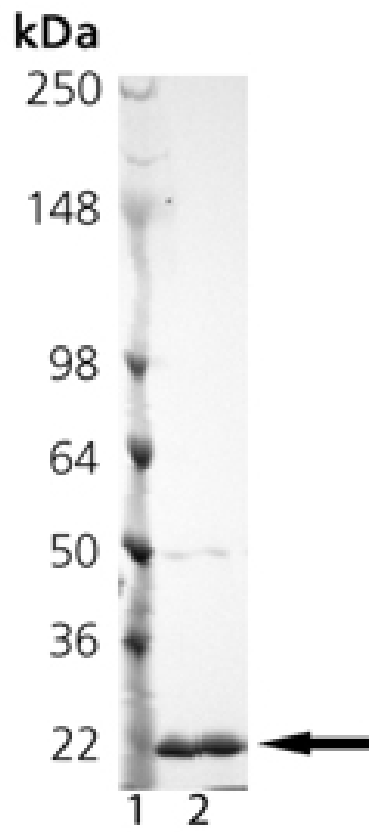
## Ordering Information

[Order Online »](#)

BML-SE722-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 2  $\mu$ g TRAIL.

## Handling & Storage

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 at 0.1 and 1.0µg/ml, in primary cultured mouse hepatocytes, in the presence of cycloheximide (5µg/ml) or actinomycin D (0.5µg/ml). Induces apoptosis in Jurkat cells (a TRAIL-sensitive human cell line) at 1µg/ml, within 3 hours. Inhibits secondary clonal expansion of mouse CD<sup>8+</sup> T cells *in vitro* (50ng/ml and greater).

**Formulation** Liquid. In PBS containing 0.1mM DTT.

**MW** 23 kDa

**Purity** ≥95% (SDS-PAGE)

**Purity Detail** Purified by multi-step chromatography.

**Source** Produced in *E. coli*. Soluble mouse TRAIL (aa 118-291).

**UniProt ID** P50591

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)