

# TRAIL, soluble (human) (recombinant) set

Citations: 20

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-850-018-KI01	1Set
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Apo-2L, TNFSF 10, CD253
Application	FUNC
Application Notes	Induces apoptosis in a concentration range of 1-100ng/ml if applied with the cross-linking enhancer.
Concentration	TRAIL: 0.1mg/ml after reconstitutionEnhancer: 1mg/ml after reconstitution
Contents	10µg of TRAIL, Soluble (human) (recombinant) (Prod. No. ALX-522-003) and 2×50µg of Enhancer for Ligands (Prod. No. ALX-804-034).
Formulation	Lyophilized. Contains PBS.
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute TRAIL with 100µl and each vial of enhancer with 50µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in <i>E. coli</i> . The extracellular domain of human TRAIL (Apo-2L) (aa 95-281) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG <sup>®</sup> -tag.
Specific Activity	ED50: 10ng/ml (Jurkat cells)
Specificity	TRAIL binds to human and mouse TRAIL receptors and human osteoprotegerin (OPG).

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)