

TRAIL (soluble) (human), (recombinant)

Citations: 24

[View Online »](#)

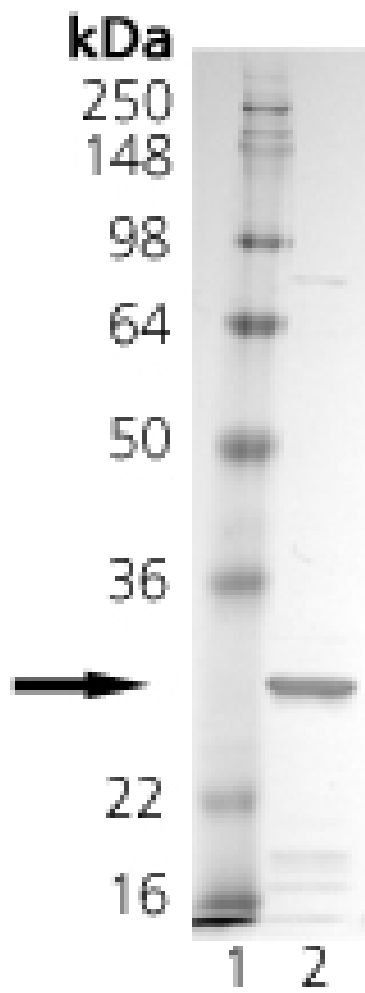
Ordering Information

[Order Online »](#)

ALX-522-003-C010	10µg
ALX-522-003-3010	SuperPack - 3x10µg

Manuals, SDS & CofA

[View Online »](#)



SDS PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g TRAIL.

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 Notes	ELISA: binds to human TRAIL receptors and human osteoprotegerin (OPG) at 10-100 ng/ml.
Biological Activity	Induces apoptosis if applied with the cross-linking enhancer (Prod. No. ALX-804-034); see also TRAIL, Soluble (human) (recombinant) Set (Prod. No. ALX-850-018). Note: rhsTRAIL does not significantly induce apoptosis in the absence of an enhancer.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~28kDa
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl 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 (aa 95-281) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.

Specificity

Binds to human and mouse TRAIL receptors and human osteoprotegerin (OPG).

Technical Info / Product Notes

Historical data has shown that Fas induces apoptosis in a concentration range of 1-100ng/ml if applied with the cross-linking enhancer (Prod. No. ALX-804-034).

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)