

TWEAK (soluble) (human), (recombinant)

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-021-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

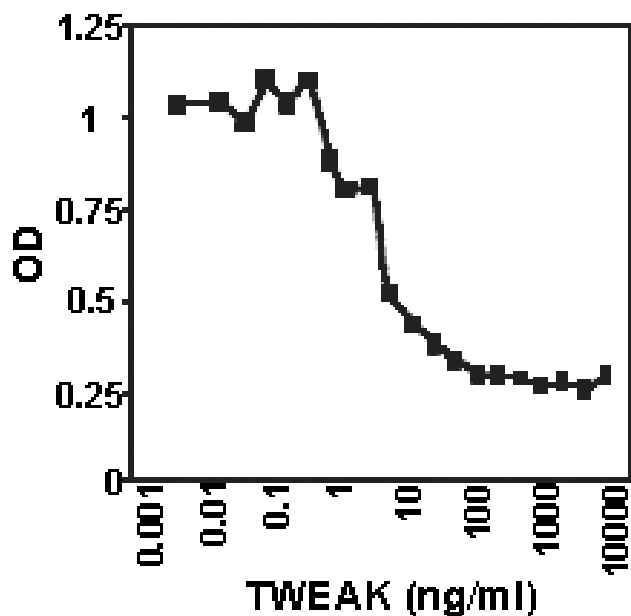


Figure: Apoptosis of TWEAK sensitive cells.

Recombinant, human TWEAK exerts its full biological activity in a concentration range of 50-500ng/ml on Kym-1 cells.**Method:** Kym-1 cells (20'000 cells in 50µl RPMI medium containing 10% fetal calf serum) were incubated with the indicated concentrations of rhsTWEAK for 24 hours at 37°C in a 96 well plate. Concentrations of rhsTWEAK required to kill target cells may vary depending on cell type studied. Cell viability was determined using a PMS/MTS-based cell proliferation assay kit.

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-3L, TNFSF 12
Application Notes	ELISA: binds to FN14 receptor at 10-100 ng/ml.
Biological Activity	Kills TWEAK sensitive cells.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~18-20kDa (SDS-PAGE).
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 TWEAK (aa 106-249) is fused at the N-terminus to a linker peptide (11 aa) and a FLAG [®] -tag.
Specificity	Binds to human and mouse Fn14 (TWEAK-R).

Technical Info / Product Notes

Historical data has shown that TWEAK has a pro-apoptotic effect on several target cell lines, and it induces angiogenesis and proliferation in endothelial cells.

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

O43508

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)