

KillerTRAIL™ Protein (soluble) (human), (recombinant) (BULK)

Citations: 8

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
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 a concentration range of 10-100ng/ml. Note: This protein does not require a cross-linking enhancer for its potent biological activity. For cell lines that require extensive cross-linking of the TRAIL-Rs for killing (e.g. Jurkat) use SUPERKILLERTRAIL [®] (Prod. No. ALX-201-115).
Endotoxin Content	<10EU/μg protein (LAL test; Bio Whittaker).
Formulation	Liquid. In 20mM HEPES, pH 7.4, containing 150mM sodium chloride, 0.005% Tween 20, 1% sucrose, 1mM DTT.
Purity	≥98% (SDS-PAGE)
Purity Detail	Affinity purified.
Reconstitution	For a special storage and dilution buffer see KillerTRAIL™ Storage and Dilution Buffer (Prod. No. ALX-505-005).
Source	Produced in <i>E. coli</i> . The extracellular domain of human TRAIL (aa 95-281) is fused at the N-terminus to a His-tag and a linker peptide.
Specific Activity	ED ₅₀ : 20ng/ml (BJAB cells)
Specificity	Binds to human and mouse TRAIL receptors and osteoprotegerin (OPG).

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)