

TRAIL (human) monoclonal antibody (2E5)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-2960	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

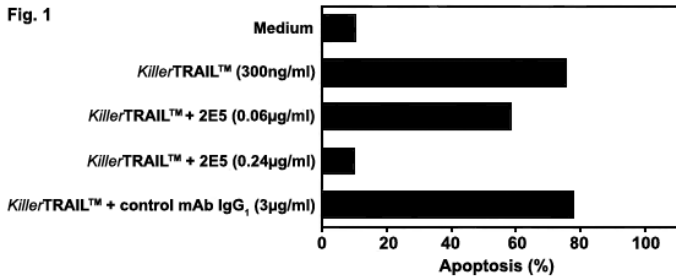
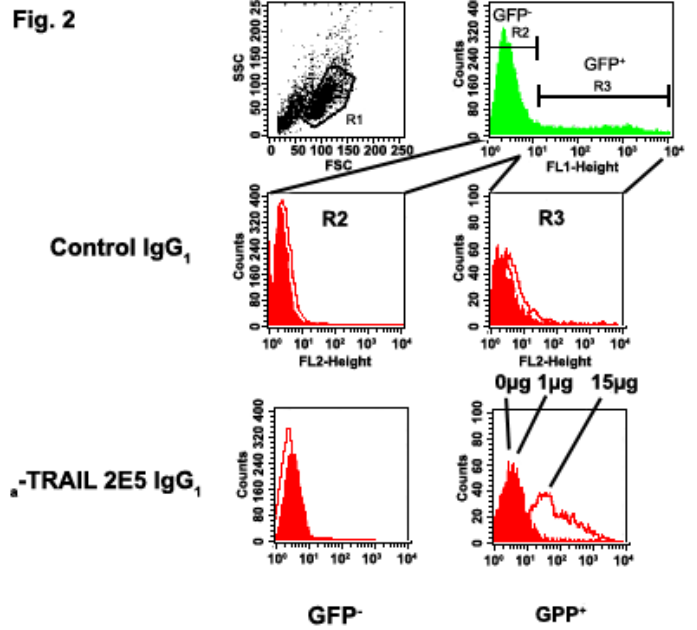


Figure 1: Apoptosis induced in Jurkat T leukemia cells by *KillerTRAIL*TM (Prod. No. [ALX-201-073](#)) is completely blocked by mouse MAb to TRAIL (human) (2E5) (Prod. No. ALX-804-296) in a concentration range of 0.5-1.5µg/ml. Concentration of MAb 2E5 required for inhibition may vary depending of the cell type studied and the concentration of soluble recombinant human TRAIL used. The neutralizing activity of MAb 2E5 has been confirmed with various sources of soluble recombinant human TRAIL including *KillerTRAIL*TM. **Method:** Jurkat cells (5×10^4) were treated with *KillerTRAIL*TM (300ng/ml) in the presence or absence of the MAb to TRAIL (human) (2E5) or a control IgG1 mouse monoclonal antibody. After 16 hours apoptosis was evaluated by forward/sideward scatter (FSC/SSC) combined with propidium iodide (PI) staining analyzed by flow cytometry.



TRAIL-207 + M2. Effect of mAbs

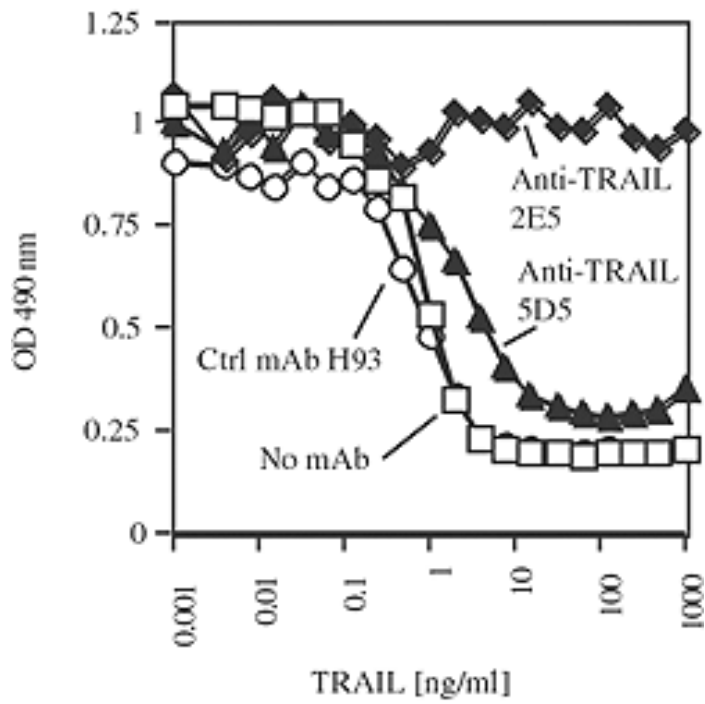


Figure: Neutralizing effect of MAb to TRAIL (human) (2E5) on TRAIL-induced apoptosis (TRAIL, Soluble (human) (recombinant) Set (Prod. No. ALX-850-018); containing Enhancer for Ligands).

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Apo-2L, TNFSF 10, CD253

Application Flow Cytometry, FUNC

Application Notes Functional Application (excellent): neutralizes human TRAIL mediated killing in a concentration range of 0.5-1.5µg/ml.
Only detects the native protein. Does not work in WB, ICC, IHC or other applications where fixation is necessary.

Clone 2E5

Crossreactivity Does not cross-react with mouse TRAIL.

Formulation Liquid. In PBS, pH 7.4. Contains no preservatives.

GenBank ID U57059

Gene/Protein Identifier 8743 (Entrez GeneID)

Host Mouse

Immunogen Recombinant human soluble TRAIL (aa 95-281).

Isotype IgG1

Purity Detail Protein A-affinity purified. Sterile filtered.

Species Reactivity Human

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)