## TRAIL (human) monoclonal antibody (2E5)

This antibody is covered by our Worry-Free Guarantee.

Citations: 11

View Online »

**Ordering Information** 

Order Online »

ALX-804-296-C100

100µg

Manuals, SDS & CofA

**View Online »** 

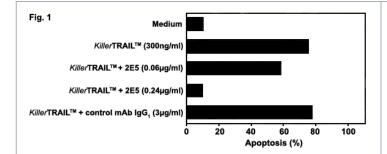
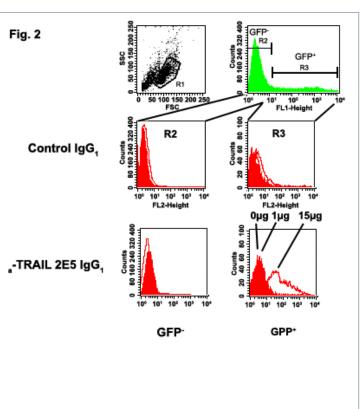
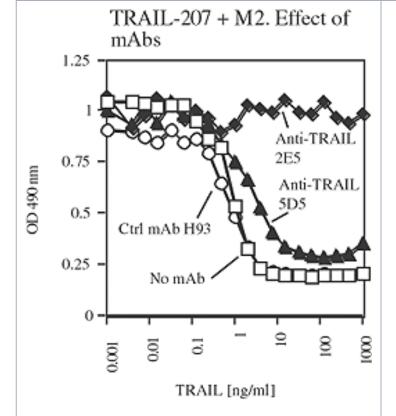


Figure 1: Apoptosis induced in Jurkat T leukemia cells by KillerTRAIL<sup>TM</sup> (Prod. No. ALX-201-073) is completely blocked by mouse MAb to TRAIL (human) (2E5) (Prod. No. ALX-804-296) in a concentration range of 0.5-1.5µg/ml. Concentration of MAb 2E5 required for inhibition may vary depending of 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<sup>TM</sup>. **Method:** Jurkat cells (5×10<sup>4</sup>) were treated with *Killer* TRAIL<sup>TM</sup> (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.""""





**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.

inc or other applications where fixation is necess

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).

lsotype lgG1

Purity Detail Protein A-affinity purified. Sterile filtered.

Species Reactivity Human

UniProt ID P50591

Last modified: May 29, 2024



info-

eu@enzolifesciences.com