

## TRAIL

Recombinant human protein expressed in E. coli cells

**Catalog # T880-40N**

Lot # 1123-6

### Product Description

Human active form TRAIL (114-281) was expressed in E. coli cells and purified by proprietary chromatographic techniques. The protein accession number is [NP\\_003801.1](#).

### Gene Aliases

TL2; APO2L; CD253; Apo-2L; TNFSF10

### Endotoxin Level

Less than 0.1 ng/µg (IEU/µg) of TRAIL.

### Formulation

Recombinant protein stored in 1 X PBS, 10% glycerol.

### Storage and Stability

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

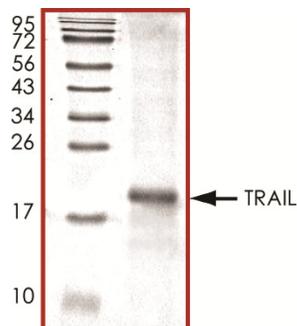
### Scientific Background

Trail or TNF-related apoptosis inducing ligand is a cytokine that belongs to the tumor necrosis factor (TNF) ligand family which function as prominent mediators of immune regulation and the inflammatory response. TRAIL induces significant cell death in human colon polyps and providing a potentially selective approach for colorectal cancer chemoprevention by targeting APC-deficient cells for apoptosis (1). TRAIL plays a crucial role in hepatic cell death and hepatic inflammation (2).

### References

- Zhang, L.. et.al: Chemoprevention of colorectal cancer by targeting APC-deficient cells for apoptosis. Nature 464: 1058-1061, 2010.
- Zheng, S.-J. et.al: Critical roles of TRAIL in hepatic cell death and hepatic inflammation. J. Clin. Invest. 113: 58-64, 2004.

### Purity



The purity was determined to be **>95%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel. Approx. MW **20.5 kDa (19523.91 Daltons)** by MW calculation.

## TRAIL

Recombinant human protein expressed in E. coli cells

Catalog Number T880-40N

Specific Lot Number 1123-6

Purity	>95%
Concentration	0.2 µg/µl
Stability	2yr at -20°C from date of shipment
Storage & Shipping	Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**