

## Anti-Trail

Rabbit Polyclonal Antibody

**Catalog # T880-63R**

Lot # O2109-3

### Cited Applications

WB, ELISA, IHC

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the TRAIL protein

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

TRAIL antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human TRAIL.

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

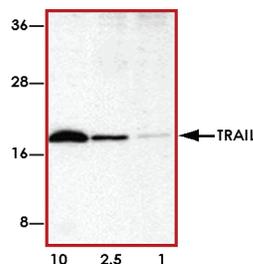
### Scientific Background

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).

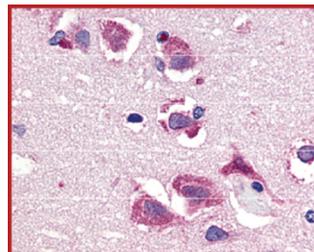
### References

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

### Sample Data



Western blot analysis of Trail in HeLa cell lysate containing 10, 2.5, or 1 ng of recombinant protein containing extracellular domain of Trail with Trail antibody at 1 µg/ml.



Immunohistochemistry of Trail in human brain tissue with Trail antibody at 20 µg/ml.

## Anti-Trail

Rabbit Polyclonal Antibody

Catalog Number

T880-63R

Specific Lot Number

O2109-3

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.**