

## TRAF1 Protein

Full-length recombinant protein expressed in Sf9 cells

**Catalog # T56-30G**

Lot # G1333-3

### Product Description

Recombinant full-length human TRAF1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_005658](#).

### Gene Aliases

EBI6, MGC:10353

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal 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.

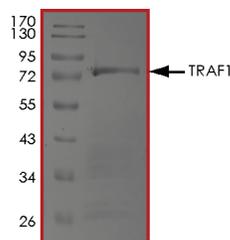
### Scientific Background

Tumor necrosis factor receptor-associated factor 1 (TRAF1) is a member of the TNF receptor-associated factor (TRAF) family. Unlike other members of the TRAF family, it lacks the N-terminal RING/zinc-finger domain. As an adapter protein, TRAF1 regulates the activation of NF-kappa-B and JNK and plays a role in the regulation of cell survival and apoptosis. TRAF1 has been shown as a substrate of PKN1 kinase activity (1). TRAF1 inhibits TNFR2-induced proteasomal degradation of TRAF2 and relieves TNFR1-induced activation of NF- $\kappa$ B from the inhibitory effect of TNFR2 (2).

### References

1. Kato T. Jr. et al: Negative regulation of constitutive NF-kappaB and JNK signaling by PKN1-mediated phosphorylation of TRAF1. *Genes Cells* 13:509-520(2008)
2. Wicovsky A. et al: Tumor necrosis factor receptor-associated factor-1 enhances proinflammatory TNF receptor-2 signaling and modifies TNFR1-TNFR2 cooperation. *Oncogene* 28:1769-1781(2009)

### Purity



The purity of TRAF1 was determined to be **>85%** by densitometry. Approx. MW **74kDa**.

## TRAF1 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog #	T56-30G
Specific Lot #	G1333-3
Purity	>85%
Concentration	0.1 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal 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.**