

## FADD Protein

Full-length recombinant human protein expressed in E. coli cells

**Catalog # F02-30H**

Lot # F553-3

### Product Description

Recombinant full-length human FADD was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM\\_003824](#).

### Gene Aliases

GIG3, MORT1, MGC8528

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

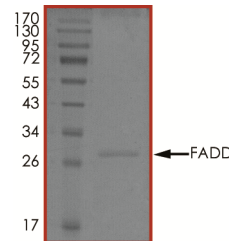
### Scientific Background

FADD or Fas-Associated protein with Death Domain is an adaptor molecule that mediates death signaling by the Fas-receptor, tumor necrosis factor receptor and TRAIL-receptor. FADD binds to these receptors via the C-terminus Death Domain which then unmask the N-terminal effector domain of FADD thereby allowing it to recruit caspase-8 and activate the cysteine protease cascade leading to *apoptosis* (1). Cells lacking FADD are defective in intracellular double-stranded RNA (dsRNA)-activated gene expression, including production of type I (alpha/beta) interferons and are thus very susceptible to viral infection (2).

### References

1. Sheikh, M.S et al: *The FADD is going nuclear*. *Cell Cycle*, 2004; 2 (4): 346-7.
2. Balachandran, S. et al: *A FADD-dependent innate immune mechanism in mammalian cells*. *Nature* 432: 401-405, 2004.

### Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **27kDa**.

## FADD Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog #	F02-30H
Lot #	F553-3
Purity	>95%
Concentration	0.2 µg/µl
Stability	1 yr 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.**