

BID Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # B59-30G

Lot # V155-3

Product Description

Recombinant full-length human BID was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC036364](#).

Gene Aliases

FP497, MGC42355, MGC15319

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°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.

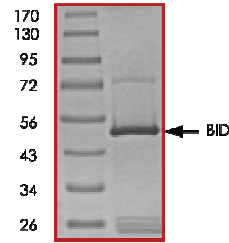
Scientific Background

BID is a BH3 interacting death domain that heterodimerizes with either agonist BAX or antagonist BCL2 (1). BID is a member of the BCL-2 family of cell death regulators and is a mediator of mitochondrial damage induced by caspase-8 (CASP8). BID initiates apoptosis by binding to regulatory sites on prosurvival BCL2 proteins to directly neutralize their function. Multiple alternatively spliced transcript variants of BID have been found, but the full-length nature of some variants has not been defined. BID together with Cathepsins play an important role in the actions of Camptothecin on breast cancer cells (2).

References

- Hayakawa, A. et al; Bid truncation mediated by caspases-3 and -9 in vinorelbine-induced apoptosis. *Apoptosis*. 2008;13(4):523-30.
- Lamparska-Przybylska, M. et al; Cathepsins and BID are involved in the molecular switch between apoptosis and autophagy in breast cancer MCF-7 cells exposed to camptothecin. *J Physiol Pharmacol*. 2005 Jun;56 Suppl 3:159-79.

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **52kDa**.

BID Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	B59-30G
Specific Lot Number	V155-3
Purity	>85%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°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
www.signalchem.com

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