

Anti-BID

Rabbit Polyclonal Antibody

Catalog # B59-363BR

Lot # O2109-30

Cited Applications

WB, ELISA

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

Specificity

Recognizes the BID protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BID antibody was raised against a peptide corresponding to 14 amino acids near the carboxy terminus of human BID

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

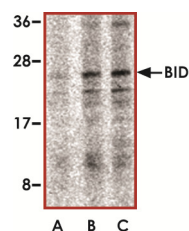
Scientific Background

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

1. Hayakawa, A. et al.; Bid truncation mediated by caspases-3 and -9 in vinorelbine-induced apoptosis. *Apoptosis*. 2008;13(4):523-30.
2. Lamparska-Przybysz, 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 PhysiolPharmacol*. 2005 Jun;56Suppl 3:159-79.

Sample Data



Western blot analysis of BID in mouse lung cell lysates with BID antibody at (A) 0.5, (B) 1, and (C) 2 ug/ml.

Anti-BID

Rabbit Polyclonal Antibody

Catalog Number

B59-363BR

Specific Lot Number

O2109-30

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
www.signalchem.com

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