

Anti-BAD

Rabbit Polyclonal Antibody

Catalog # B03-363R

Lot # O2109-27

Cited Applications

WB, ELISA, IF, IHC

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

Specificity

Recognizes the BAD protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BAD antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human BAD

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

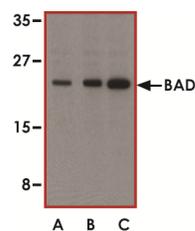
Scientific Background

BAD is a member of the BCL-2 family of proteins that are known to be regulators of programmed cell death. BAD is a pro-apoptotic protein that forms a heterodimer complex with BCL-xL and BCL-2 which reverses the prosurvival activity of these proteins (1). The proapoptotic activity of BAD is regulated through its phosphorylation and this inhibits the pro-apoptosis function of BAD. Protein kinases such as AKT, RAF and RSK1 can phosphorylate BAD and RSK1-induced phosphorylation of BAD at ser112 suppresses BAD-mediated apoptosis in neurons. BAD inhibits G(1) to S phase transition in MCF7 breast cancer cells and overexpression of BAD inhibits cell growth as well as cyclin D1 expression (2).

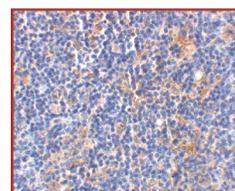
References

1. Yang E, et al: Bad, a heterodimeric partner for Bcl-X(L) and Bcl-2, displaces Bax and promotes cell death. Cell 80: 285-291, 1995.
2. Fernando R, et al: Breast cancer cell proliferation is inhibited by BAD: regulation of cyclin D1. J Biol Chem. 2007 Sep 28;282(39):28864-73.

Sample Data



Western blot analysis of BAD in T24 cell lysates with BAD antibody at (A) 0.5, (B) 1, and (C) 2 µg/ml.



Immunohistochemical staining of rat thymus using BAD at 2 µg/ml.

Anti-BAD

Rabbit Polyclonal Antibody

Catalog Number

B03-363R

Specific Lot Number

O2109-27

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.