

Anti-BAG1

Rabbit Polyclonal Antibody

Catalog # B05-363BR

Lot # O2109-54

Cited Applications

WB, ELISA, IF, IHC

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

Specificity

Recognizes the BAG1 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

BAG1 antibody was raised against a 14 amino acid synthetic peptide from near the amino terminus of human BAG1

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

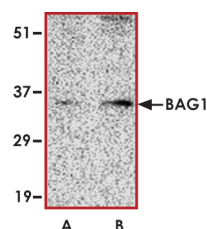
Scientific Background

BAG1 (also known as BCL2-associated athanogene) is a membrane protein rich in glutamic acid residues that binds to BCL2 and blocks apoptosis or programmed cell death (1). The BAG1-BCL2 complex enhances the anti-apoptotic effects of BCL2 and represents a link between growth factor receptors and anti-apoptotic mechanisms (2). Overexpression of BAG1 in 3T3 fibroblasts prevents apoptosis in the presence of low serum. BAG1 has also been shown to interact with activated glucocorticoid, androgen, estrogen and progesterone receptors. Binding to these receptors by BAG1 is dependent on receptor activation.

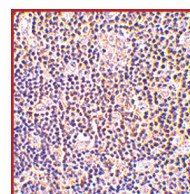
References

1. Tang, S.C. BAG-1, an anti-apoptotic tumour marker.". IUBMB Life, 2003; 53 (2): 99-105.
2. Clemo, N.K. et al: The role of the retinoblastoma protein (Rb) in the nuclear localization of BAG-1: implications for colorectal tumour cell survival.". Biochem. Soc. Trans.2005; 33 (Pt 4): 676-8.

Sample Data



Western blot analysis of BAG1 in PC-3 cell lysate with BAG1 antibody at (A) 1 and (B) 2 µg/ml.



Immunohistochemistry of BAG1 in human lymph node tissue with BAG1 antibody at 2 µg/ml

Anti-BAG1

Rabbit Polyclonal Antibody

Catalog Number

B05-363BR

Specific Lot Number

O2109-54

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.