

BAG1 Protein

Recombinant human protein expressed in *E. coli* cells

Catalog # **B05-31G**

Lot # D122-4

Product Description

Recombinant human BAG1 (72-end) was expressed in *E. coli* cells using an N-terminal GST tag. The gene accession number is [NM_004323](#).

Gene Aliases

RAP46

Formulation

Recombinant protein stored in 50mM MOPS, pH 7.5, 150mM NaCl, 150mM Imidazole, 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

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.

Purity



BAG1 Protein

Recombinant human protein expressed in *E. coli* cells

Catalog Number	B05-31G
Specific Lot Number	D122-4
Purity	>85%
Concentration	0.2 $\mu\text{g}/\mu\text{l}$
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.