

BAD (Ab-134) Antibody [#B0822]

Catalog Number:	B0822
Concentration:	1mg/ml
Swiss-Prot No.:	Q92934
Other Names:	BAD, BBC6, BCL2L8, Bcl- XL/Bcl-2 associated death promoter, Bcl-2 binding component 6, Bcl-2-like 8 protein, Bcl-xL/Bcl-2 associated death promoter, Bcl2-antagonist of cell death
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human BAD around the phosphorylation site of serine 134 (P-K-S ^P -A-G).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BAD (Ab-134) antibody detects endogenous levels of total BAD protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Ottillie S., J. Biol. Chem. 272:30866-30872(1997). Wang H.-G., Cell 87:629-638(1996). Petros A.M., Protein Sci. 9:2528-2534(2000).