

## BAD Antibody [#B0821]

<b>Catalog Number:</b>	B0821
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q92934
<b>Other Names:</b>	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
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from Internal of human BAD.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	BAD antibody detects endogenous levels of total BAD protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Wang H.-G., Cell 87:629-638(1996). Ottilie S., J. Biol. Chem. 272:30866-30872(1997). The MGC Project Team, Genome Res. 14:2121-2127(2004).