



BAD (Phospho-Ser112) Antibody

Catalog Number 11067

(For Research Use Only)

- Catalog Number:** 11067-1, 11067-2, 11067
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Storage/Stability: Store at -20°C/1 year
Immunogen: The antiserum was produced against synthesized phosphopeptide derived from human BAD around the phosphorylation site of serine 112 (H-S-S^P-Y-P).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity/Sensitivity: BAD (phospho-Ser112) antibody detects endogenous levels of BAD only when phosphorylated at serine 112.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:4000
Swiss-Prot No. : Q92934
References: Zhang B, et al. (2004). Mol Cell Biol.24 (14): 6205-6214.
Rice PL, et al. (2003). Cancer Res.63 (3): 616-620.
Wang XQ, et al. (2001). J Biol Chem.276 (48): 44504-44511.

