

BAD (Ab-155) Antibody [#B7019]

- Catalog Number:** B7019-1, B7019-2, B7019
- 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:** This antiserum was produced against synthesized non-phosphopeptide derived from human BAD around the phosphorylation site of serine 155.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** BAD (Ab-155) antibody detects endogenous levels of total BAD protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:10000
- Swiss-Prot No. :** Q92934
- References:** Moon EY, et al. (2003). Blood.101 (10): 4122-4130.