

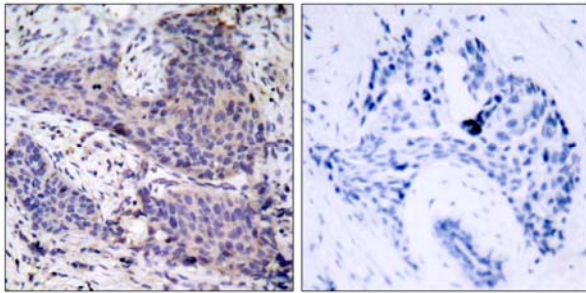


BAD (Ab-136) Antibody

Catalog Number 21063

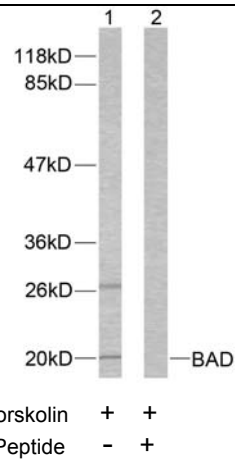
(For Research Use Only)

- Catalog Number:** 21063-1, 21063-2, 21063
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 serine136.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: BAD (Ab-136) antibody detects endogenous levels of total BAD protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
Swiss-Prot No. : Q92934
References: Wang XQ, et al. (2001). J Biol Chem.276 (48): 44504-44511.
Lee YI, et al. (2001). J Biol Chem.276 (20): 16969-16977.
Maiti D, et al. (2001). J Biol Chem.276 (1): 329-333.



Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using BAD (Ab-136) antibody (#21063).



Western blot analysis of extracts from 293 cells using BAD (Ab-136) antibody (#21063).