

## BID (Ab-78) Antibody [#B0461]

<b>Catalog Number:</b>	B0461
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P55957
<b>Other Names:</b>	BH3 interacting domain death agonist
<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 non-phosphopeptide derived from human BID around the phosphorylation site of serine 78 (S-E-S <sup>P</sup> -Q-E).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	BID (Ab-78) antibody detects endogenous levels of total BID protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Wang K., Genes Dev. 10:2859-2869(1996). Footz T.K., Genomics 51:472-475(1998). Renshaw S.A., J. Biol. Chem. 279:2846-2855(2004).