

BIK (Ab-33) Antibody [#B0053]

- Catalog Number:** B0053-1, B0053-2, B0053
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q13323
- All Names:** Apoptosis inducer NBK, BIKLK, BIP1, BP4, Bcl-2 interacting killer, NBK
- 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 non-phosphopeptide derived from human BIK around the phosphorylation site of threonine 33 (G-M-T-D-S).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** BIK (Ab-33) antibody detects endogenous levels of total BIK protein.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
IF: 1:500~1:1000 ELISA: 1:10000
- References:** Sulekha Verma, J. Biol. Chem., Feb 2001; 276: 4671 - 4676.
Yan M. Li, Cancer Res., Nov 2003; 63: 7630 - 7633.
Malti Nikrad, Mol. Cancer Ther., Mar 2005; 4: 443 - 449.
Feng Dong, Infect. Immun., Mar 2005; 73: 1861 - 1864.
Hiroo Ueno, Mol. Biol. Cell, Feb 2000; 11: 735.
Hidenori Hase, J. Biol. Chem., Nov 2002; 277: 46950 - 46958.
Yasuyuki Hirashima, J. Biol. Chem., Jul 2003; 278: 26793 - 26802.