

## Zap-70 (Ab-493) Antibody [#B7256]

- Catalog Number:** B7256-1, B7256-2, B7256
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P43403
- All Names:** 70 kDa zeta-associated protein, SRK, Syk-related tyrosine kinase, ZA70, ZAP-70, ZAP70, kinase ZAP70
- Form of Antibody:** 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human Zap-70 around the phosphorylation site of tyrosine 493 (S-Y-Y<sup>P</sup>-T-A).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Zap-70 (Ab-493) antibody detects endogenous levels of total Zap-70 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** IHC: 1:50~1:100 ELISA: 1:20000
- References:** Chan AC, et al. (1995) EMBO J; 14(11): 2499-2508.  
Kong G, et al. (1996) Mol Cell Biol; 16(9): 5026-5035.  
Salomon AR, et al. (2003) Proc Natl Acad Sci USA; 100(2): 443-448.