

Zap-70 (Ab-319) Antibody [#B7255]

- Catalog Number:** B7255-1, B7255-2, B7255
- 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 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 Zap-70 around the phosphorylation site of tyrosine 319 (S-P-Y^P-S-D).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Zap-70 (Ab-319) antibody detects endogenous levels of total Zap-70 protein. This antibody may cross-react with Syk when phosphorylated at equivalent sites.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
- References:** Salomon AR, et al. (2003) Proc Natl Acad Sci USA; 100(2): 443-448.
Ku GM, et al. (2001) EMBO J; 20(3): 457-465.
Tang J, et al. (1999) Proc Natl Acad Sci U S A; 96(17): 9775-9780.
Zhao Q, et al. (1996) Mol Cell Biol; 16(12): 6765-6774.
Williams BL, et al. (1999) EMBO J; 18(7): 1832-1844.
Adjali O, et al. (2005) J Clin Invest; 115(8): 2287-2295.