



ASK1(Phospho-Ser966) Antibody

Catalog Number 11179

(For Research Use Only)

- Catalog Number:** 11179-1, 11179-2, 11179
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: The antiserum was produced against synthesized phosphopeptide derived from human ASK1 around the phosphorylation site of Serine 966 (S-I-S^P-L-P).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity/Sensitivity: ASK1(Phospho-Ser966) Antibody detects endogenous levels of ASK1 only when phosphorylated at Serine 966.
Reactivity: Human, Mouse
Applications: IHC: 1:50~1:100 ELISA: 1:20000
Swiss-Prot No. : Q99683
References: Zhang W, et al. (2005) J Biol Chem. 280(19): 19036-19044.
Fujii K, et al. (2004) Oncogene. 23(29):5099-5104.
Goldman EH, et al. (2004) J Biol Chem 2004 Mar 12; 279(11): 10442-10449.
Zhang L, et al. (1999) Proc Natl Acad Sci U S A. 96(15): 8511-8515.

