



ASK1(Phospho-Ser83) Antibody

Catalog Number 11178

(For Research Use Only)

- Catalog Number:** 11178-1, 11178-2, 11178
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 83.
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-Ser83) Antibody detects endogenous levels of ASK1 only when phosphorylated at Serine 83.
Reactivity: Human
Applications: WB: 1:500~1:1000 ELISA: 1:20000
Swiss-Prot No. : Q99683
References: Mabuchi S, et al. (2004) *Endocrinology*. 145(1): 49-58.
Yuan ZQ, et al. (2003) *J Biol Chem*. 278(26): 23432-23440.
Kim AH, et al. (2001) *Mol Cell Biol*. 21(3): 893-901.

