



cdc25C (Phospho-Ser216) Antibody

Catalog Number 11118

(For Research Use Only)

- Catalog Number:** 11118-1, 11118-2, 11118
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : P30307
All Names: CDC25M1, Dual specificity phosphatase Cdc25C, M-phase inducer phosphatase 3, MPIP3
All Sites: Human: Ser216
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 cdc25C around the phosphorylation site of serine 216 (S-P-S^P-M-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: cdc25C (phospho-Ser216) antibody detects endogenous levels of cdc25C only when phosphorylated at serine 216.
Reactivity: Human
Applications: IHC: 1:50~1:100 ELISA: 1:10000
References: Toyoshima-Morimoto F. et al. (2002) EMBO Rep. 3(4): 341-348.
Ferguson AM. et al. (2005) Mol Cell Biol. 25(7): 2853-2860.
Donzelli M. et al. (2003) EMBO Rep. 4(7): 671-677.
Chen F. et al. (2002) Proc Natl Acad Sci U S A. 99(4): 1990-1995.



