



CDC2 (Phospho-Tyr15) Antibody

Catalog Number 11244

(For Research Use Only)

- Catalog Number:** 11244-1, 11244-2, 11244
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : P06493
All Names: p34 protein kinase; Cyclin-dependent kinase 1; CDK1
All Sites: Human: Tyr15; Mouse: Tyr15; Rat:Tyr15
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 CDC2 around the phosphorylation site of tyrosine 15 (G-T-Y^P-G-V).
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: CDC2 (Phospho-Tyr15) antibody detects endogenous levels of CDC2 only when phosphorylated at tyrosine 15.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:10000
References: Jeremy P. H. Chow, et al. (2003) J. Biol. Chem; 278: 40815-40828.
NJ Wells, et al. (1999) J. Cell Sci; 112: 3361.
P Jin, et al. (1996) J. Cell Biol; 134: 963.
KM Pickham and DJ Donoghue (1994) Mol. Biol. Cell; 5: 587.

