



CDC2 (Ab-15) Antibody

Catalog Number 21236

(For Research Use Only)

Catalog Number: 21236-1, 21236-2, 21236

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No. : P06493

All Names: p34 protein kinase; Cyclin-dependent kinase 1; CDK1

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 immunogen.

Specificity/Sensitivity CDC2 (Ab-15) antibody detects endogenous levels of total CDC2 protein.

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.

