

RCC1 Antibody [#C0314]

Catalog Number: C0314-1, C0314-2, C0314

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

Swiss-Prot No. : P18754

All Names: CHC1, Cell cycle regulatory protein, Regulator of chromosome condensation

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 peptide derived from human RCC1.

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity RCC1 antibody detects endogenous levels of total RCC1 protein.

Reactivity: Human

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:5000

References: M Dasso, Mol. Cell. Biol., Aug 1992; 12: 3337 - 3345.
Hoi Yeung Li, J. Cell Biol., Mar 2003; 160: 635.
Michael E. Nemerlut, Science, May 2001; 292: 1540.
Ian Cushman, Mol. Biol. Cell, Jan 2004; 15: 245.