



## Cortactin (Ab-421) Antibody

Catalog Number 21263

(For Research Use Only)

- Catalog Number:** 21263-1, 21263-2, 21263  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Swiss-Prot No. :** Q14247  
**All Names:** Amplexin, CTTN, EMS1, Oncogene EMS1, SRC8, Src substrate cortactin  
**Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), 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 non-phosphopeptide derived from human Cortactin around the phosphorylation site of tyrosine 421 (P-I-Y<sup>P</sup>-E-D).  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** Cortactin (Ab-421) antibody detects endogenous levels of total Cortactin protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000 ELISA: 1:20000  
**References:** Du, Y. et al. (1998) Mol. Cell. Biol. 18, 5838-5851.  
Naisbitt, S. et al. (1999) Neuron 23, 569-582.  
Rodrigo, J. P. et al. (2000) Clin. Cancer Res. 6, 3177-3182.  
Weed, S.A. and Parsons, J.T. (2001) Oncogene 20, 6418-6434.

