



## cdc25C (Ab-216) Antibody

Catalog Number 21145

(For Research Use Only)

- Catalog Number:** 21145-1, 21145, 21145  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Swiss-Prot No. :** P30307  
**All Names:** CDC25M1, Dual specificity phosphatase Cdc25C, M-phase inducer phosphatase 3, MPIP3  
**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 cdc25C around the phosphorylation site of serine 216 (S-P-S<sup>P</sup>-M-P).  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** cdc25C (Ab-216) antibody detects endogenous levels of total cdc25C protein.  
**Reactivity:** Human  
**Applications:** IHC: 1:50~1:100 ELISA: 1:20000  
**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.



