



## CaMKII (Ab-286) Antibody

Catalog Number 21279

(For Research Use Only)

- Catalog Number:** 21279-1, 21279-2, 21279  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Swiss-Prot No. :** Q9UQM7  
**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 CaMKII around the phosphorylation site of threonine 286 (Q-E-T<sup>P</sup>-V-D).  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** CaMKII (Ab-286) antibody detects endogenous levels of total CaMKII protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000  
**References:** Pak JH, et al. Proc Natl Acad Sci U S A. 2000 Oct 10; 97(21): 11232-11237  
Hudmon A, et al. J Cell Biol. Author manuscript; available in PMC 2006 May 7  
Miller P, et al. PLoS Biol. 2005 Apr; 3(4): e107  
Runyan JD, et al. Learn Mem. 2005 Mar; 12(2): 103-110.

