



CaMKII (Phospho-Thr286) Antibody

Catalog Number 11287

(For Research Use Only)

- Catalog Number:** 11287-1, 11287-2, 11287
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : Q9UQM7
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 CaMKII around the phosphorylation site of threonine 286 (Q-E-T^P-V-D).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity/Sensitivity: CaMKII (phospho-Thr286) antibody detects endogenous levels of CaMKII only when phosphorylated at threonine 286.
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.

