

Product Information



CaMKII Monoclonal Antibody (Clone 6G9)

Item No. 10011437

Contents:	This vial contains protein G-purified mouse IgG at a concentration of 1 mg/ml in PBS, pH. 7.4, containing 50% glycerol and 0.09% sodium azide.
Synonym:	Calcium/Calmodulin-dependent Protein Kinase II
Antigen:	Rat recombinant CaMKII
Host:	Mouse, clone 6G9
Isotype:	IgG ₁
Cross Reactivity:	(+) Mouse, rat, and bovine CaMKII. Detects ~50-60 kDa protein, corresponding to the apparent molecular mass of CaMKII α -subunit on SDS-PAGE immunoblots. Both the phosphorylated and non-phosphorylated forms are recognized.
Stability:	≥ 1 year at -20°C
Applications:	Immunofluorescence, immunoprecipitation, western blot (WB), immunohistochemistry, and ELISA. The recommended starting concentration for WB is 0.1 μ g/ml; 0.1 μ g/ml of CaMKII antibody was needed for detection of CaMKII in 20 μ g of rat brain tissue. Optimal working dilutions for other applications should be determined empirically.

CaMKII is an important member of the calcium/calmodulin-activated protein kinase family, functioning in neural synaptic stimulation and T-cell receptor signaling.^{1,2} CaMKII is expressed in many different tissues but is specifically found in neurons of the forebrain and its mRNA is found within the dendrites and the soma of the neuron. The CaMKII that is found in the neurons consist of two subunits of 52 (termed α genes) and 60 kDa (β genes). CaMKII has catalytic and regulatory domains, as well as an ATP-binding domain, and a consensus phosphorylation site.³⁻⁷ The binding of Ca²⁺/calmodulin to its regulatory domain releases its auto inhibitory effect and activates the kinase.⁸ This kinase activation results in autophosphorylation at threonine 286 (Thr²⁸⁶).⁸

The threonine phosphorylation state of CaMKII can be regulated through protein phosphatase 1 (PP1)/protein kinase A (PKA). Whereas PP1 dephosphorylates phospho-CaMKII at Thr²⁸⁶, PKA prevents this dephosphorylation.⁹ Autophosphorylation also enables CaMKII to attain an enhanced affinity for NMDA receptors in postsynaptic densities.¹⁰⁻¹²

References

1. Hughes, K., Edin, S., Antonsson, Å., *et al. J. Biol. Chem.* **276(38)**, 36008-36013 (2001).
2. Barria, A., Muller, D., Derkach, V., *et al. Science* **276**, 2042-2045 (1997).
3. Bennett, M.K. and Kennedy, M.B. *Proc. Natl. Acad. Sci. USA* **84**, 1794-1798 (1987).
4. Brocke, L., Srinivassan, M., and Schulman, H. *J. Neurosci.* **15(10)**, 6797-6808 (1995).
5. Nghiem, P., Saati, S.M., Martens, C.L., *et al. J. Biol. Chem.* **268(8)**, 5471-5479 (1983).
6. Edman, C.F. and Schulman, H. *Biochim. Biophys. Acta* **1221(1)**, 89-101 (1994).
7. Tombes, R.M. and Krystal, G.W. *Biochim. Biophys. Acta* **1355(3)**, 281-292 (1997).
8. Means, A.R. *Mol. Endocrinol.* **14(1)**, 4-13 (2000).
9. Makhinson, M., Chotiner, J.K., Watson, J.B., *et al. J. Neurochem.* **19(7)**, 2500-2510 (1999).
10. Strack, S. and Colbran, R.J. *J. Biol. Chem.* **273(33)**, 20689-20692 (1998).
11. Leonard, A.S., Lim, I.A., Hemsworth, D.E., *et al. Proc. Natl. Acad. Sci. USA* **96**, 3239-3244 (1999).
12. Shen, K. and Meyer, T. *Science* **284**, 162-167 (1999).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10011437

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 12/19/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com