

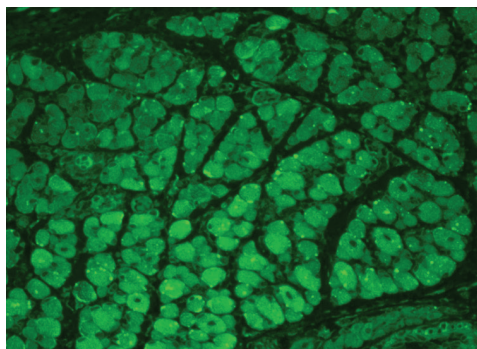
# PRODUCT INFORMATION

## CaMKII (Phospho-Thr<sup>286</sup>/Thr<sup>287</sup>) Monoclonal Antibody (Clone 22B1) Item No. 10011438

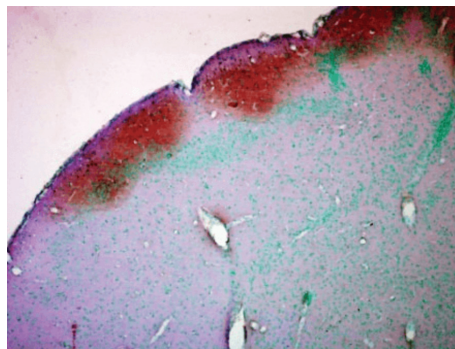
### Overview and Properties

<b>Contents:</b>	This vial contains 25 or 100 µg of protein G-affinity purified monoclonal antibody.
<b>Synonym:</b>	Calcium/Calmodulin-dependent Protein Kinase II
<b>Immunogen:</b>	Synthetic peptide
<b>Cross Reactivity:</b>	(+) Human, mouse, rat; other species not tested
<b>Uniprot No.:</b>	P11275
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, pH. 7.4, with 50% glycerol and 0.09% sodium azide
<b>Concentration:</b>	1 mg/ml
<b>Clone:</b>	22B1
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG1
<b>Applications:</b>	ELISA, Immunocytochemistry (ICC), Immunofluorescence (IF), Immunohistochemistry (IHC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:1,000 for ICC, IF, and WB, and 1:100 for IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Immunohistochemical labeling of mouse backskin tissue. Tissue was fixed with Bouin's Fixative and embedded in paraffin and incubated with CaMKII (Phospho-Thr<sup>286</sup>/Thr<sup>287</sup>) Monoclonal Antibody (Clone 22B1) at a 1:100 dilution for one hour at room temperature. Then tissues were labeled with FITC Goat Anti-Mouse (green) at 1:50 for one hour at room temperature. Positive staining was localized to muscle, hair follicle, epidermis.



Immunohistochemical staining of formalin-fixed human colon carcinoma tissue using CaMKII (Phospho-Thr<sup>286</sup>/Thr<sup>287</sup>) Monoclonal Antibody (Clone 22B1) at a 1:5,000 dilution for 12 hours at 4°C. Following incubation with Biotin Goat Anti-Mouse at a dilution of 1:2,000 for one hour at room temperature and counterstained with 200 µl Mayer Hematoxylin (purple/blue) nuclear stain for two minutes at room temperature.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/20/2023

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

CaMKII is an important member of the calcium/calmodulin-activated protein kinase family, functioning in neural synaptic stimulation and T-cell receptor signaling.<sup>1,2</sup> CaMKII is expressed in many different tissues but is specifically found in the 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 consists of two subunits of 52 (termed  $\alpha$  genes) and 60 kDa ( $\beta$  genes). CaMKII has catalytic and regulatory domains, as well as an ATP-binding domain, and a consensus phosphorylation site.<sup>3-7</sup> The binding of  $\text{Ca}^{2+}$ /calmodulin to its regulatory domain releases its autoinhibitory effect and activates the kinase.<sup>8</sup> This kinase activation results in autophosphorylation at threonine 286 (Thr<sup>286</sup>).<sup>8</sup> The threonine phosphorylation state of CaMKII can be regulated through protein phosphatase 1 (PP1)/protein kinase A (PKA). Whereas PP1 dephosphorylates phospho-CaMKII at Thr<sup>286</sup>, PKA prevents this dephosphorylation.<sup>9</sup> Autophosphorylation also enables CaMKII to attain an enhanced affinity for NMDA receptors in postsynaptic densities.<sup>10-12</sup>

## References

1. Hughes, K., Edin, S., Antonsson, Å., *et al.* Calmodulin-dependent kinase II mediates T cell receptor/CD3- and phorbol ester-induced activation of I $\kappa$ B kinase. *J. Biol. Chem.* **276**(38), 36008-36013 (2001).
2. Barria, A., Muller, D., Derkach, V., *et al.* Regulatory phosphorylation of AMPA-type glutamate receptors by CaM-KII during long-term potentiation. *Science* **276**(7), 2042-2045 (1997).
3. Bennett, M.K. and Kennedy, M.B. Deduced primary structure of the  $\beta$  subunit of brain type II  $\text{Ca}^{2+}$ /calmodulin-dependent protein kinase determined by molecular cloning. *Proc. Natl. Acad. Sci. USA* **84**(7), 1794-1798 (1987).
4. Brocke, L., Srinivasan, M., and Schulman, H. Development and regional expression of multifunctional  $\text{Ca}^{2+}$ /calmodulin-dependent protein kinase isoforms in rat brain. *J. Neurosci.* **15**(10), 6797-6808 (1995).
5. Nghiem, P., Saati, S.M., Martens, C.L., *et al.* Cloning and analysis of two new isoforms of multifunctional  $\text{Ca}^{2+}$ /calmodulin-dependent protein kinase. Expression in multiple human tissues. *J. Biol. Chem.* **268**(8), 5471-5479 (1983).
6. Edman, C.F. and Schulman, H. Identification and characterization of  $\delta_B$ -CaM kinase and  $\delta_C$ -CaM kinase from rat heart, two new multifunctional  $\text{Ca}^{2+}$ /calmodulin-dependent protein kinase isoforms. *Biochim. Biophys. Acta* **1221**(1), 89-101 (1994).
7. Tombes, R.M. and Krystal, G.W. Identification of novel human tumor cell-specific CaMK-II variants. *Biochim. Biophys. Acta* **1355**(3), 281-292 (1997).
8. Means, A.R. Regulatory cascades involving calmodulin-dependent protein kinases. *Mol. Endocrinol.* **14**(1), 4-13 (2000).
9. Makhinson, M., Chotiner, J.K., Watson, J.B., *et al.* Adenylyl cyclase activation modulates activity-dependent changes in synaptic strength and  $\text{Ca}^{2+}$ /calmodulin-dependent kinase II autophosphorylation. *J. Neurochem.* **19**(7), 2500-2510 (1999).
10. Strack, S. and Colbran, R.J. Autophosphorylation-dependent targeting of calcium/calmodulin-dependent protein kinase II by the NR2B subunit of the N-methyl-D-aspartate receptor. *J. Biol. Chem.* **273**(33), 20689-20692 (1998).
11. Leonard, A.S., Lim, I.A., Hemsworth, D.E., *et al.* Calcium/calmodulin-dependent protein kinase II is associated with the N-methyl-D-aspartate receptor. *Proc. Natl. Acad. Sci. USA* **96**(6), 3239-3244 (1999).
12. Shen, K. and Meyer, T. Dynamic control of CaMKII translocation and localization in hippocampal neurons by NMDA receptor stimulation. *Science* **284**(5411), 162-166 (1999).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM