

CaMKII polyclonal antibody

CaMKII (calmodulin-dependent kinase II) is an enzyme which is activated by increases in intracellular Ca^{2+} ion concentration and it has been proposed to be pivotal in regulating synaptic strength and maturation of synapses during development. This process is thought to be critical in memory and learning and in establishing the specificity of synaptic connections. There are over two dozen alternative splice variants of CaMKII which are encoded by four genes, α , β , γ , and δ with apparent molecular masses of 50-60 kDa. CaMKII is widely distributed in many tissues, but is highly expressed in brain. Several studies demonstrate that CaMKII is a multifunctional enzyme which modulates the synaptic strength by binding to a subunit of NMDA receptors and promoting the phosphorylation of this NMDA receptor subunit. CaMKII also regulates DLG localization at synapses by co-localization with DLG in the same protein complex. Experimental data suggest that CaMKII is critically involved in the development of morphine tolerance as well as dependence and inhibition of this enzyme may have some therapeutic benefit in the treatment of opiate tolerance and dependence. It also has been demonstrated that δ CaMKII isozyme is down-regulated in human tumor cells indicating a role for δ CaMKII isozymes in cellular differentiation. Changes in δ CaMKII isozyme expression pattern in human hearts during heart failure suggest that CaMKII is important for regulation of heart function. This immunoaffinity purified antibody detects proteins of 50-60 kDa, corresponding to apparent molecular mass of CaMKII isoforms on SDS-PAGE immunoblots, in samples from human, mouse, rat, bovine, hamster, guinea pig, chicken and rabbit origins. Recombinant rat calmodulin-dependent protein kinase II α subunits expressed in sf9 insect cells are also detected. As the sequences of the rat α , δ , and γ isoforms are conserved over this amino terminus region, this antibody is expected to recognize the δ , γ and α isoforms. This antibody is not expected to cross-react with the β isoform. Proteins of unknown identity, ~40, 45 and 90 kDa, may be detected on immunoblot analysis with some lysates.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

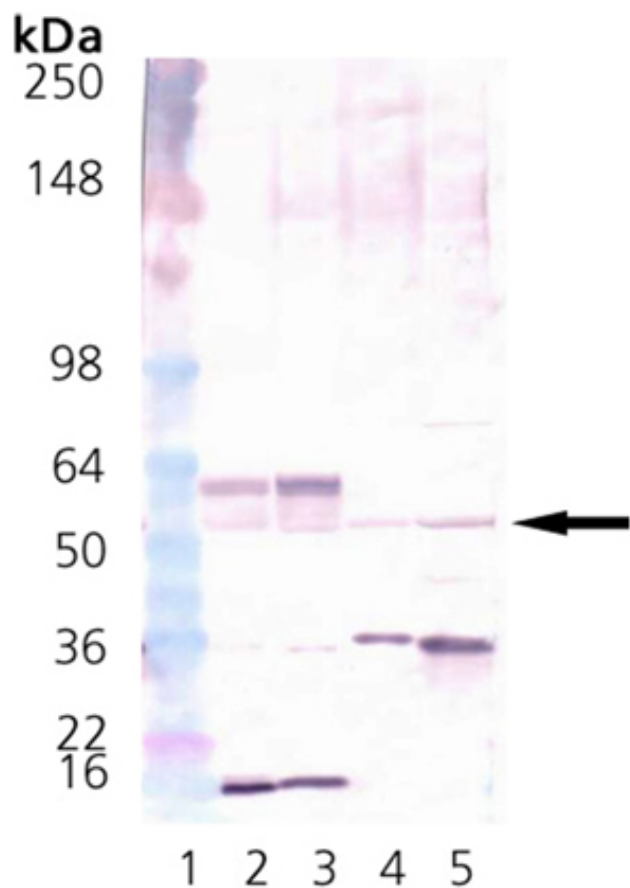
[Order Online »](#)

ADI-KAP-CA002-D	50µg
-----------------	------

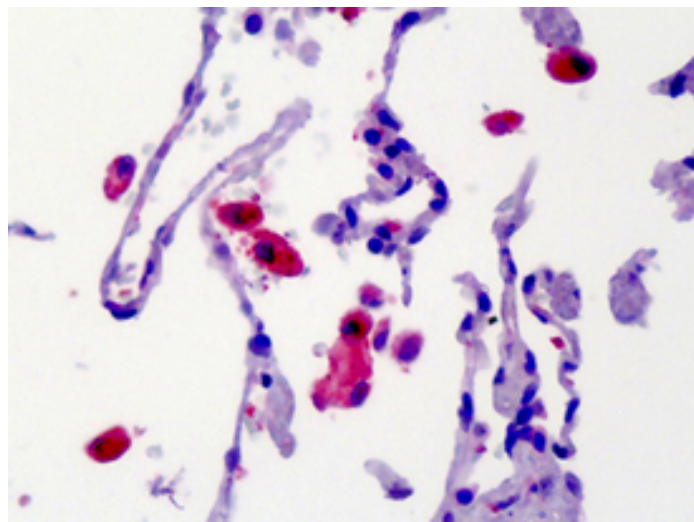
ADI-KAP-CA002-F	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of CaMKII: Lane 1: MWM, Lane 2: Brain (mouse), (tissue extract) (Prod No. ADI-LYTMB100) Lane 3: Brain (rat), (tissue extract) (Prod No. ADI-LYT-RB100) Lane 4: HeLa, (cell lysate) (Prod No. ADI-LYC-HL100) Lane 5: Hs-67 cell lysate.



IHC Analysis: Human lung tissue stained with CaMKII, pAb at 10µg/ml.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Calcium/calmodulin-dependent protein kinase II

Application IHC (PS), WB

Application Notes Detects a band of ~50-60kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

GenBank ID J05072

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of rat CaMKII δ .

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (Colorimetric, 1:500) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Chicken, Guinea pig, Hamster, Human, Mouse, Rabbit, Rat

UniProt ID P15791

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)