CaMKII polyclonal antibody

CaMKII (calmodulin-dependent kinase II) is an enzyme which is activated by increases in intracellular Ca²⁺ 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, a, b, g, and d 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 d CaMKII isozyme is down-regulated in human tumor cells indicating a role for d CaMKII isozymes in cellular differentiation. Changes in d 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 a subunits expressed in sf9 insect cells are also detected. As the sequences of the rat a, d, and g isoforms are conserved over this amino terminus region, this antibody is expected to recognize the d, g and a isoforms. This antibody is not expected to crossreact with the b 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 »

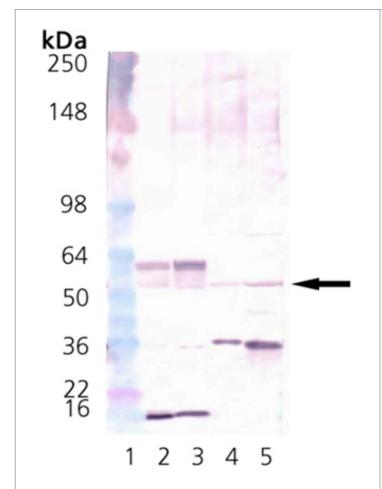
ADI-KAP-CA002-D

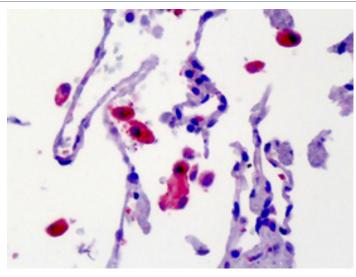
50µg

ADI-KAP-CA002-F 200μg

Manuals, SDS & CofA

View Online »





IHC Analysis: Human lung tissue stained with CaMKII, pAb at $10\mu g/ml$.

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.

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com