

CAMKII Monoclonal Antibody [6G9]

(Cat. No. A-0456)

Background

CaM-kinase II (CAMK2) is a member of the NMDAR signaling complex in excitatory synapses and may regulate NMDAR-dependent potentiation of the AMPAR and synaptic plasticity. Moreover, it is a prominent kinase in the central nervous system that may function in long-term potentiation and neurotransmitter release.

Description

Mouse monoclonal antibody to CAMKII, Clone 6G9

Formulation

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

Immunogen

Rat CaMKII (partially purified).

Specificity

Mouse, Rat

Isotype

IgG1

Purification

Protein G- affinity purified.

Storage

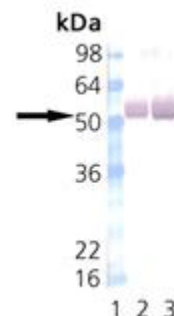
Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

Calcium/calmodulin-dependent protein kinase II

Application

IHC, WB (Colorimetric, 1:1000)



▲ **Fig. 1.** Western Blot Analysis of Calmodulin-Dependent Protein Kinase II: Lane 1: MWM, Lane 2: Mouse Brain, Lane 3: Rat Brain.

Ordering Information

Products

CAMKII Monoclonal Antibody [6G9]

Size

50 µg

Cat. No.

A-0456-050

This product is for research purposes only. Not intended for use in diagnostic procedures.