# PKA (CT) polyclonal antibody

PKA (Protein Kinase A), also known as cAMP-dependent protein kinase, phosphorylates serine or threonine residues in target protein in response to elevated levels of cAMP. Inactive PKA exists as a tetrameric protein composed of two regulatory (R) subunits and two catalytic (C) subunits. Activation occurs when two cAMP molecules bind to each R subunit which causes a conformational change that releases the active C subunits.

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

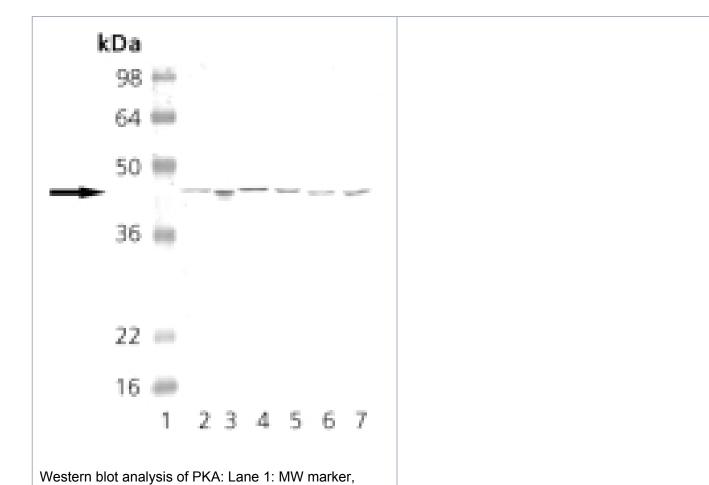
## **Ordering Information**

**Order Online** »

ADI-KAS-PK017-D	50µg
ADI-KAS-PK017-F	200μg

Manuals, SDS & CofA

**View Online »** 



Lane 2: CHO-K1, Lane 3: Mouse Brain, Lane 4: Rat Brain, Lane 5: HeLa, Lane 6: 3T3, Lane 7: PC-12.

#### **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 Protein Kinase A

**Application** WB

**Application Notes**Detects a band of ~40kDa by Western blot.

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

azide.

**Host** Rabbit

Immunogen Synthetic peptide corresponding to a portion of human

PKA.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, colorimetric)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, Dog, Guinea pig, Hamster, Human, Monkey,

Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

**UniProt ID** P17612 (Cα subunit), P22694 (Cβ subunit), P22162 (Cγ

subunit), P10644 (RIα subunit), P31321 (RIβ subunit),

P13861 (RIIa subunit), P31323 (RIIB subunit)

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com