

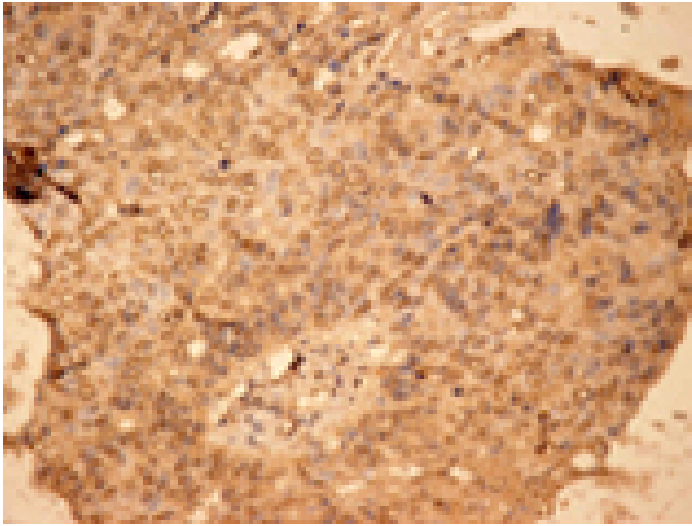
# PKC polyclonal antibody

Protein Kinase C (PKC) is a large superfamily of serine/threonine kinases that mediate essential cellular signals required for activation, proliferation, differentiation and survival. There are at least ten PKC isotypes that are closely related in structure but that have distinct patterns of tissue distribution and function. The PKC isotypes can be subdivided into three classes based on primary structure and biochemical properties. These are: classical or conventional PKC isotypes (cPKC), novel PKC isotypes (nPKC) and atypical PKC isotypes (aPKC).

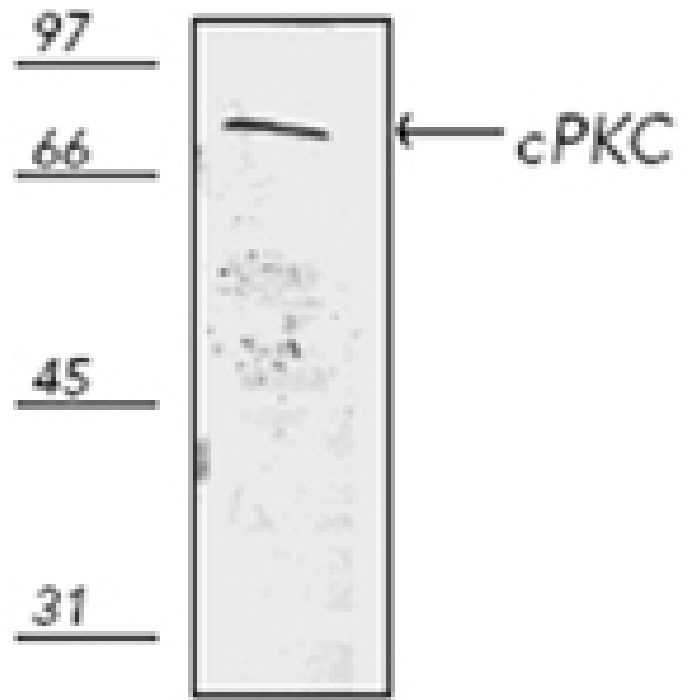
This antibody is covered by our [Worry-Free Guarantee](#).

**Manuals, SDS & CofA**

[View Online »](#)



Immunohistochemistry analysis of human lung cancer tissue with PKC pAb.



Western blot analysis of human thymus HS67 cell lysate, probed with PKC pAb.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	protein kinase C
Application	IHC, IP, WB
Application Notes	Detects a band of ~67-83kDa by Western blot.
Formulation	Liquid. In borate buffered saline, pH 8.2.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human PKC. The sequence is completely conserved between the $\alpha$ , $\beta$ , and $\delta$ isoforms of PKC.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50)Immunoprecipitation (12.5 $\mu$ g/ml)Western Blot (2 $\mu$ g/ml, ECL)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, Dog, Guinea pig, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep
UniProt ID	P17252 (PKCalpha)

### 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)

Last modified: May 29, 2024