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.

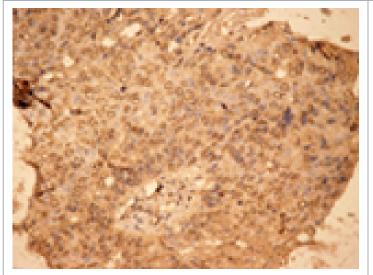
Ordering Information

Order Online »

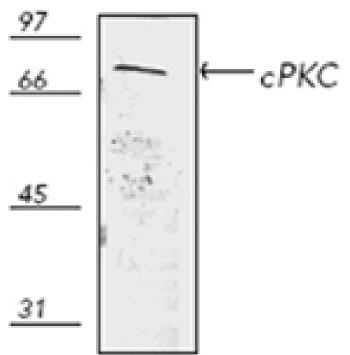
ADI-KAP-PK003-E 100μg

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

 α , β , and δ isoforms of PKC.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:50)Immunoprecipitation

(12.5µg/ml)Western Blot (2µg/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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com