

Anti-PKAc

Rabbit Polyclonal Antibody

Catalog # P51-63R

Lot # D316-5

Cited Applications

- Western blot (1:250)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the PKAc α , PKAc β and PKAc γ protein

Cross Reactivity

- Western blot human PKAc α , PKAc β and PKAc γ

PKAc α , PKAc β and PKAc γ from other species may also be detectable

Host

Rabbit

Immunogen

Protein Code: P17612

Formulation

TBS pH 7.2, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

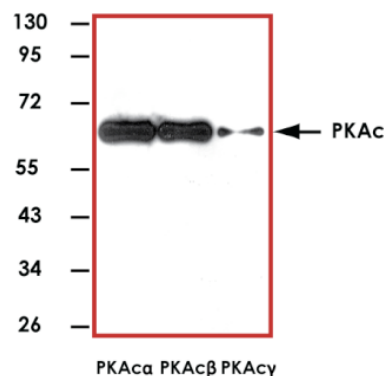
The second messenger cyclic AMP (cAMP) activates cAMP-dependent protein kinase (PKA or cAPK) in mammalian cells and controls many cellular mechanisms such as gene transcription, ion transport and protein phosphorylation (1). Inactive PKA is a heterotetramer composed of a regulatory subunit (R) dimer and a catalytic subunit (C) dimer, and in this inactive state, the pseudosubstrate sequences on the R subunits block the active sites on the C subunits. Three C subunit isoforms (- α -, - β - and - γ -) and two families of regulatory subunits (RI and RII) with distinct cAMP binding properties have been identified. Upon binding of cAMP to the R subunits, the autoinhibitory contact is eased and active monomeric C subunits are released.

References

1. Montminy M: Transcriptional regulation by cyclic AMP. *Annu Rev Biochem*, 1997;66:807-22.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-(604)-232-4601 or by email: orders@signalchem.com
www.signalchem.com

Sample Data



Representative western blot with Anti-PKAc (1:250) using 20ng of PKAc α , PKAc β and PKAc γ recombinant protein.

Anti-PKAc

Rabbit Polyclonal Antibody

Catalog Number

P51-63R

Specific Lot Number

D316-5

Purification

Affinity purified using Protein G

Concentration

0.25 μ g/ μ L

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.