

PRKAR1A Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # P57-30H

Lot # M370-1

Product Description

Recombinant full-length human PRKAR1A Protein was expressed in Sf9 insect cells using a C-terminal His tag. The gene accession number is [NM_002734](#).

Gene Aliases

CAR; CNC; CNC1; DKFZp779L0468; MGC17251; PKR1; PPNAD1; PRKAR1; TSE1

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

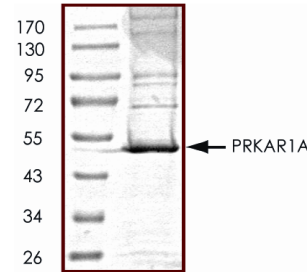
Scientific Background

PRKAR1A or the cAMP-dependent protein kinase regulatory subunit type I alpha, is part of the type I PKA holoenzyme. PRKAR1A is found to be a tissue-specific extinguisher that down-regulates the expression of seven liver genes in hepatoma x fibroblast hybrids (1). Mutations in PRKAR1A gene cause Carney complex (CNC) and PRKAR1A can fuse to the RET protooncogene by gene rearrangement and form the thyroid tumor-specific chimeric oncogene known as PTC2 (2). A nonconventional nuclear localization sequence (NLS) has been found for this protein which suggests a role in DNA replication via the protein serving as a nuclear transport protein for the second subunit of the Replication Factor C (RFC40).

References

1. Bossis I, et al: PRKAR1A: normal and abnormal functions. *Endocrinology* 145: 5452-5458, 2004.
2. Casey M, et al: Mutations in the protein kinase A R1-alpha regulatory subunit cause familial cardiac myxomas and Carney complex. *J. Clin. Invest.* 106: R31-R38, 2000.

Purity



The purity of PRKAR1A was determined to be **>85%** by densitometry. Approx. MW **51kDa**.

PRKAR1A Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number P57-30H

Specific Lot Number M370-1

Purity	>85%
Concentration	0.2 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.