

PRKY Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P82-35G

Lot # T1039-2

Product Description

Recombinant human PRKY (40-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC074852](#).

Gene Aliases

(None)

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 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.

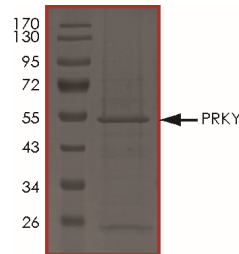
Scientific Background

PRKY or protein kinase, Y-linked, pseudogene is a novel serine/threonine subtype of protein kinase that appears to encode a protein related to the catalytic subunit of the cAMP-dependent protein kinases which are key players in the cellular responses to the second messenger cAMP. PRKY gene is classified as a transcribed pseudogene because it has lost a coding exon that results in all transcripts being candidates for nonsense-mediated decay (NMD) and unlikely to express a protein (1). Abnormal XY interchange between a novel isolated protein kinase gene, PRKY, and its homologue, PRKX, accounts for one third of all (Y+)XX males and (Y-)XY females (2).

References

- Schiebel, K.et.al: FISH localization of the human Y-homolog of protein kinase PRKX (PRKY) to Yp11.2 and two pseudogenes to 15q26 and Xq12-q13. Cytogenet. Cell Genet. 76: 49-52, 1997.
- Schiebel, K.et.al: Abnormal XY interchange between a novel isolated protein kinase gene, PRKY, and its homologue, PRKX, accounts for one third of all (Y+)XX males and (Y-)XY females. Hum. Molec. Genet. 6: 1985-1989, 1997.

Purity



The purity of PRKY protein was determined to be **>75%** by densitometry. Approx. MW **53 kDa**.

PRKY Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	P82-35G
Lot #	T1039-2
Purity	>75%
Concentration	0.05 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.