

TP53RK Protein

Recombinant full-length protein expressed in Sf9 cells

Catalog # T29-34G

Lot # J677-3

Product Description

Recombinant full-length human TP53RK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The TP53RK gene accession number is [NM_033550](#).

Gene Aliases

BUD32; C20orf64; dJ101A2; Nori-2; Nori-2p

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM 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.

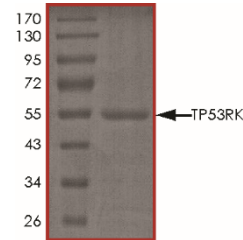
Scientific Background

TP53RK or TP53 regulating kinase and also called PRPK (p53-related protein kinase) is the human homologue of yeast Bud32. TP53RK belongs to a small subfamily of atypical protein kinases. TP53RK is expressed in interleukin-2-activated cytotoxic T-cells, epithelial tumor cell lines, and the testes and lower expression of TP53RK is found in heart, kidney, and spleen (1). PRPK interacts with CGI-121 and co-precipitation of p53 with PRPK is inhibited by adding recombinant CGI-121 in vitro. CGI-121 may act as an inhibitor of the PRPK-p53 interaction (2). TP53RK is activated by another kinase, Akt/PKB, which phosphorylates TP53RK at Ser250.

References

1. Abe, Y.et.al: Cloning and characterization of a p53-related protein kinase expressed in interleukin-2-activated cytotoxic T-cells, epithelial tumor cell lines, and the testes. J. Biol. Chem. 276: 44003-44011, 2001.
2. Miyoshi, A.et.al: Identification of CGI-121, a novel PRPK (p53-related protein kinase)-binding protein. Biochem. Biophys. Res. Commun. 303: 399-405, 2003.

Purity



The purity of TP53RK was determined to be **>90%** by densitometry, approx. MW **55 kDa**.

TP53RK Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	T29-34G
Lot #	J677-3
Purity	>90%
Concentration	0.1 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.