

PTK9 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P97-34G

Lot # X466-6

Product Description

Recombinant human PTK9 (2-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The PTK9 gene accession number is [NM_002822](#).

Gene Aliases

Twinfilin 1; TWF1; A6

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.

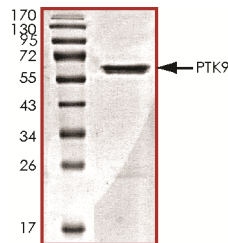
Scientific Background

PTK9 is an actin monomer-binding protein conserved from yeast to mammals which has been localized to cortical G-actin-rich structures that may be regulated by the small GTPase RAC1. PTK9 has tyrosine kinase activity similar to that of recombinant FGFR2, with optimal activity over pH 6.5 to 7.4 and a preference for manganese over magnesium as a divalent cofactor (1). PTK9 plays a central role in regulation of actin dynamics and cell motility in lymphoma cell homeostasis in vivo and represent as a bona fide lymphoma drug target (2).

References

1. Beeler. et.al: Prokaryotic expression cloning of a novel human tyrosine kinase. *Molec. Cell. Biol.* 14: 982-988, 1994.
2. Meacham. et.al: In vivo RNAi screening identifies regulators of actin dynamics as key determinants of lymphoma progression. *Nature Genet.* 41: 1133-1137, 2009.

Purity



The purity of PTK9 was determined to be **>95%** by densitometry, approx. MW **64 kDa**.

PTK9 Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	P97-34G
Specific Lot Number	X466-6
Purity	>95%
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.