

PTK9L (A6R) Protein

Full-length recombinant human protein expressed in sf9 cells

Catalog # P98-30G

Lot # U1775-6

Product Description

Full-length recombinant human PTK9L (A6R) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC003161](#).

Gene Aliases

TWF2, A6r, A6RP, MSTP011

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.

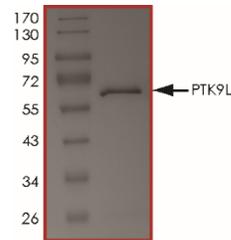
Scientific Background

Protein tyrosine kinase 9-like (A6-related protein) belongs to the actin-binding proteins ADF family, the Twinfilin subfamily. It was initially cloned by interaction using the catalytic domain of protein kinase Czeta (PKCzeta) as bait. It contains actin-binding sites and ATP-binding sites. Involved in motile and morphological processes by regulating actin polymerization.

References

1. Rohwer A, et al: Cloning, expression and characterization of an A6-related protein. Eur J Biochem. 263(2):518-25. 1999.
2. <http://www.uniprot.org/uniprot/Q6IBS0>

Purity



The purity of PTK9L (A6R) protein was determined to be **>95%** by densitometry, PTK9L (A6R) approx. MW **65kDa**.

PTK9L (A6R) Protein

Full-length recombinant human protein expressed in sf9 cells

Catalog #	P98-30G
Lot #	U1775-6
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.