

Recombinant PTK6 protein

Catalog No: 81339, 81639

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.4 µg/µl

Source: Human

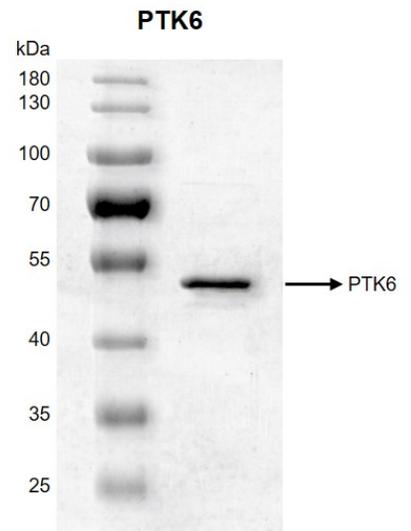
Buffer Contents: Recombinant PTK6 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: PTK6 is a non-receptor tyrosine-protein kinase which implicated in the regulation of a variety of signaling pathways and control the differentiation and maintenance of normal epithelia, as well as tumor growth. Its function seems to be context dependent and differ depending on cell type, as well as its intracellular localization. A number of potential nuclear and cytoplasmic substrates have been identified. These include the RNA-binding proteins: KHDRBS1/SAM68, KHDRBS2/SLM1, KHDRBS3/SLM2 and SFPQ/PSF; transcription factors: STAT3 and STAT5A/B and a variety of signaling molecules: ARHGAP35/p190RhoGAP, PXN/paxillin, BTK/ATK, STAP2/BKS. It also associates with a variety of proteins that are likely upstream of PTK6 in various signaling pathways, or for which PTK6 may play an adapter-like role. These proteins include ADAM15, EGFR, ERBB2, ERBB3 and IRS4. Association with EGFR, ERBB2, ERBB3 may contribute to mammary tumor development and growth through enhancement of EGF-induced signaling via BTK/AKT and PI3 kinase. PTK6 also contributes to migration and proliferation by contributing to EGF-mediated phosphorylation of ARHGAP35/ p190RhoGAP, which promotes association with RASA1/p120RasGAP, inactivating RhoA while activating RAS.

Protein Details: Recombinant PTK6 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_005966.1) with a N-terminal FLAG Tag. The molecular weight of the protein is 53.1 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant PTK6 protein gel
10% SDS-PAGE with Coomassie blue staining

MW: 53 kDa

Purity: >95%