

PTK7 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # P95-31G

Lot # J428-1

Product Description

Recombinant human PTK7 (726-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_002821](#).

Gene Aliases

CCK-4; CCK4

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.

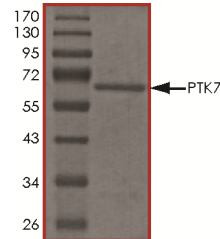
Scientific Background

PTK7 or pseudo protein tyrosine kinase 7 is a member of the receptor protein tyrosine kinase family of proteins that transduce extracellular signals across the cell membrane. PTK7 lacks detectable catalytic tyrosine kinase activity but is involved in the Wnt signaling pathway and plays a role in multiple cellular processes including polarity and adhesion (1). PTK7 interacts with β -catenin and functions upstream from glycogen synthase kinase 3. PTK7 has been isolated from colon carcinoma and was designated colon carcinoma kinase-4 (2). In conjunction with other T cell markers, PTK7 has utility as a biomarker for detecting minimal residual disease of T-ALL in the bone marrow.

References

- Lu, X.et.al: PTK7/CCK-4 is a novel regulator of planar cell polarity in vertebrates. *Nature* 430: 93-98, 2004.
- Mossie, K. et.al: Colon carcinoma kinase-4 defines a new subclass of the receptor tyrosine kinase family. *Oncogene* 11: 2179-2184, 1995.

Purity



The purity of PTK7 was determined to be **>95%** by densitometry. Approx. MW **66 kDa**.

PTK7 Protein

Recombinant human protein expressed in Sf9 cells

| | |
|---------------------|--|
| Catalog Number | P95-31G |
| Specific Lot Number | J428-1 |
| 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.