

PFTK1 (CDK14)/Cyclin Y, Active

Recombinant human proteins expressed in Sf9 cells

Catalog # P15-10G

Lot # U1535-2

Product Description

Recombinant human full-length PFTK1 (CDK14) and Cyclin Y (2-end) were co-expressed by baculovirus in Sf9 insect cells using N-terminal GST tags. The PFTK1 (CDK14) gene accession number is [NM_012395](#) and Cyclin Y is [NM_145012](#).

Gene Aliases

PFTK1: CDK14; PFTAIRE1

Cyclin Y: CCNY; C10orf9; CBCP1; CCNX; CFP1

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.

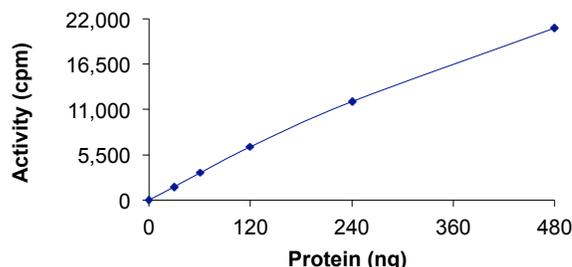
Scientific Background

PFTK1 is a member of the CDC2 - related protein kinase family which is expressed primarily in the postnatal and adult nervous system (1). PFTK1 is expressed at high levels in brain, pancreas, kidney, heart, testis, and ovary. PFTK1 amino acid residue ser119 is required for its interaction with all four 14-3-3 isoforms, while ser120 may contribute slightly to PFTK1 binding (2). PFTK1 acts as a cyclin-dependent kinase that regulates cell cycle progression and cell proliferation. Overexpression of PFTK1 predicts resistance to chemotherapy in patients with oesophageal squamous cell carcinoma.

References

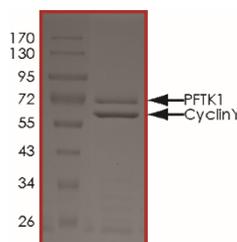
1. Yang, T. et.al: Identification and cellular localization of human PFTAIRE1. *Gene* 267: 165-172, 2001.
2. Gao, Y. et.al: A Cdc2-related protein kinase hPFTAIRE1 from human brain interacting with 14-3-3 proteins. *Cell Res.* 16: 539-547, 2006.

Specific Activity



The specific activity of PFTK1 (CDK14)/Cyclin Y was determined to be **2.7 nmol/min/mg** as per activity assay protocol.

Purity



The purity of PFTK1 (CDK14)/Cyclin Y was determined to be **>90%** by densitometry. PFTK1 (CDK14) approx. MW **72kDa** and Cyclin Y Approx. MW **64kDa**.

PFTK1 (CDK14)/Cyclin Y, Active

Recombinant human proteins expressed in Sf9 cells

Catalog #	P15-10G
Specific Activity	2.7 nmol/min/mg
Lot #	U1535-2
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.