

PCTK1 (CDK16)/CyclinY, Active

Recombinant human proteins expressed in Sf9 cells

Catalog # P09-11G

Lot # O929-2

Product Description

Recombinant human PCTK1 (CDK16) (107-end) and Cyclin Y (2-end) were co-expressed by baculovirus in Sf9 insect cells using N-terminal GST tags. The PCTK1 (CDK16) gene accession number is [NM_006201](#) and Cyclin Y is [NM_145012](#).

Gene Aliases

PCTK1 (CDK16): FLJ16665; PCTAIRE; PCTAIRE1; PCTGAIRE
CyclinY: 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.

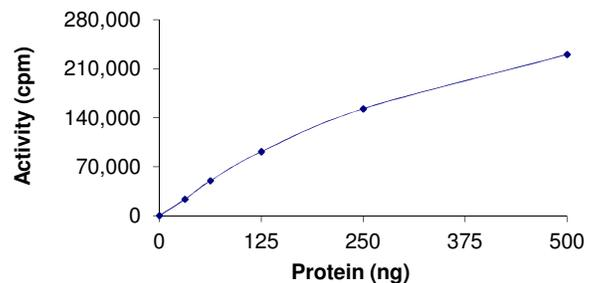
Scientific Background

PCTK1 or PCTAIRE PROTEIN KINASE 1 belongs to the cdc2/cdkx subfamily of the ser/thr family of protein kinases. PCTK1 plays a role in signal transduction cascades in terminally differentiated cells, in exocytosis, and in transport of secretory cargo from the endoplasmic reticulum (1). PCTK1 is ubiquitously expressed with the highest levels detected in the brain and testis. PCTK1 also contained the gene coding for ubiquitin-activating enzyme UBE1 that previously shown to mapped to Xp11.3 (2).

References

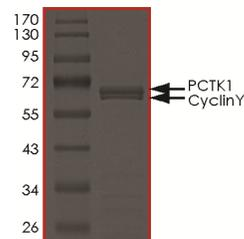
- Meyerson, M. et.al: A family of human cdc2-related protein kinases. EMBO J. 11: 2909-2917, 1992.
- Knight, J. C. et.al: Physical linkage of the cdc2-related gene (PCTK1) and the ubiquitin-activating enzyme E1 gene (UBE1) on human Xp11.3. Cytogenet. Cell Genet. 71: 155-157, 1995.

Specific Activity



The specific activity of PCTK1 (CDK16)/Cyclin Y was determined to be **40 nmol /min/mg** as per activity assay protocol.

Purity



The purity of PCTK1 (CDK16) /Cyclin Y was determined to be **>90%** by densitometry, PCTK1 (CDK16) approx. MW **66kDa** and Cyclin Y Approx. MW **64kDa**.

PCTK1 (CDK16)/CyclinY, Active

Recombinant human proteins expressed in Sf9 cells

Catalog #	P09-11G
Specific Activity	40 nmol/min/mg
Lot #	O929-2
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.

Activity Assay Protocol

Reaction Components

Active Kinase (Catalog #: P09-11G)

Active PCTK1 (CDK16)/Cyclin Y (0.1µg/µl) diluted with Kinase Dilution Buffer III (Catalog #: K23-09) and assayed as outlined in sample activity plot. (Note: these are suggested working dilutions and it is recommended that the researcher perform a serial dilution of Active PCTK1 (CDK16)/Cyclin Y for optimal results).

Kinase Dilution Buffer III (Catalog #: K23-09)

Kinase Assay Buffer I (Catalog #: K01-09) diluted at a 1:4 ratio (5X dilution) with 50ng/µl BSA solution.

Kinase Assay Buffer I (Catalog #: K01-09)

Buffer components: 25mM MOPS, pH 7. 2, 12.5mM β-glycerol-phosphate, 25mM MgCl₂, 5mM EGTA, 2mM EDTA. Add 0.25mM DTT to Kinase Assay Buffer prior to use.

[³³P]-ATP Assay Cocktail

Prepare 250µM [³³P]-ATP Assay Cocktail in a designated radioactive working area by adding the following components: 150µl of 10mM ATP Stock Solution (Catalog #: A50-09), 100µl [³³P]-ATP (1mCi/100µl), 5.75ml of Kinase Assay Buffer I (Catalog #: K01-09). Store 1ml aliquots at -20°C.

10mM ATP Stock Solution (Catalog #: A50-09)

Prepare ATP stock solution by dissolving 55mg of ATP in 10ml of Kinase Assay Buffer I (Catalog #: K01-09). Store 200µl aliquots at -20°C.

Substrate (Catalog #: M42-51N)

MBP Protein diluted in distilled H₂O to a final concentration of 1mg/ml.

Note: Histone H1 protein (Catalog #H10-54N) and Rb (C-term) protein (Catalog #R05-55G) have also been used as the substrates for this target and they showed good activity.

Assay Protocol

- Step 1.** Thaw [³³P]-ATP Assay Cocktail in shielded container in a designated radioactive working area.
- Step 2.** Thaw the Active PCTK1 (CDK16)/Cyclin Y, Kinase Assay Buffer, Substrate and Kinase Dilution Buffer on ice.
- Step 3.** In a pre-cooled microfuge tube, add the following reaction components bringing the initial reaction volume up to 20µl:
 - Component 1.** 10µl of diluted Active PCTK1 (CDK16)/Cyclin Y (Catalog #P09-11G)
 - Component 2.** 5µl of 1mg/ml stock solution of substrate (Catalog #M42-51N)
 - Component 3.** 5µl distilled H₂O (4°C)
- Step 4.** Set up the blank control as outlined in step 3, excluding the addition of the substrate. Replace the substrate with an equal volume of distilled H₂O.
- Step 5.** Initiate the reaction by the addition of 5 µl [³³P]-ATP Assay Cocktail bringing the final volume up to 25µl and incubate the mixture in a water bath at 30°C for 15 minutes.
- Step 6.** After the 15 minute incubation period, terminate the reaction by spotting 20 µl of the reaction mixture onto individual pre-cut strips of phosphocellulose P81 paper.
- Step 7.** Air dry the pre-cut P81 strip and sequentially wash in a 1% phosphoric acid solution (dilute 10ml of phosphoric acid and make a 1L solution with distilled H₂O) with constant gentle stirring. It is recommended that the strips be washed a total of 3 intervals for approximately 10 minutes each.
- Step 8.** Count the radioactivity on the P81 paper in the presence of scintillation fluid in a scintillation counter.
- Step 9.** Determine the corrected cpm by removing the blank control value (see Step 4) for each sample and calculate the kinase specific activity as outlined below.

Calculation of [³³P]-ATP Specific Activity (SA) (cpm/pmol)

Specific activity (SA) = cpm for 5 µl [³³P]-ATP / pmoles of ATP (in 5 µl of a 250 µM ATP stock solution, i.e., 1250 pmoles)

Kinase Specific Activity (SA) (pmol/min/µg or nmol/min/mg)

Corrected cpm from reaction / [(SA of ³³P-ATP in cpm/pmol)*(Reaction time in min)*(Enzyme amount in µg or mg)]*[(Reaction Volume) / (Spot Volume)]

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.