

TTK, Active

Full-length recombinant protein expressed in Sf9 cells

Catalog # T20-10G Lot # I327-3

Product Description

Recombinant full-length human TTK was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is <u>NM 003318</u>.

Gene Aliases

ESK, PYT, MPS1, MPS1L1, FLJ38280

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

TTK is a serine/threonine kinase that has been implicated in the regulation of centrosome duplication and mitotic checkpoint response. Overexpressing of dominantnegative TTK in human cell lines prevents centrosome duplication while active TTK accelerates centrosome reduplication in an osteosarcoma cell line. Disruption of TTK function leads to a combination of severe mitotic abnormalities and failure in centrosome duplication (1). TTK is required for stabilization and activation of p53 during spindle disruption. TTK phoshorylates the N-terminal domain of p53 at Thr18, and this phosphorylation disrupts the interaction with MDM2 and abrogates MDM2mediated p53 ubiquitination (2).

References

- 1. Fisk H A, et al: Human Mps1 protein kinase is required for centrosome duplication and normal mitotic progression. *Proc. Nat. Acad. Sci.* 100: 14875-14880, 2003.
- 2. Huang Y F, et al: TTK/hMps1 mediates the p53-dependent postmitotic checkpoint by phosphorylating p53 at Thr18. Mol Cell Biol. 2009 Jun;29(11):2935-44.

 Catalog #
 Aliquot Size

 T20-10G -05
 5 μg

 T20-10G -10
 10 μg

Specific Activity



The specific activity of TTK was determined to be **18 nmol** /min/mg as per activity assay protocol.

Purity



The purity of TTK was determined to be **>75%** by densitometry, approx. MW **~130kDa**.

TTK, Active

Full-length recombinant protein expressed in Sf9 cells

Catalog Number Specific Activity Specific Lot Number Purity

> Concentration Stability Storage & Shipping

T20-10G 18 nmol/min/mg I327-3 >75%

0.1 µg/µl lyr at -70°C from date of shipment 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: <u>orders@signalchem.com</u> <u>www.signalchem.com</u>

Activity Assay Protocol

Reaction Components

Active Kinase (Catalog #: T20-10G)

Active TTK $(0.1\mu g/\mu l)$ 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 TTK 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 MgC1₂, 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 200 μ l aliquots at -20°C.

Substrate (Catalog #: M42-51N)

Myelin Basic Protein (MBP) substrate diluted in distilled H_2O to a final concentration of 1mg/ml.

Assay Protocol

- Step 1. Thaw [³³P]-ATP Assay Cocktail in shielded container in a designated radioactive working area.
- Step 2. Thaw the Active TTK, 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 TTK (Catalog #T20-10G)
 - 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 [P³³]-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: <u>orders@signalchem.com</u> www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.