

Data Sheet

CDK17/Cyclin Y, GST-tags

Human, Recombinant, N-terminal GST-tags

Catalog #: 100606

Lot #: 191217

Conc.: 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 50 mM NaCl, 10 mM glutathione, 0.25 mM DTT, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Meyerson, M. *et al*, *EMBO J.* **11**: 2909-2917, 1992.
2. Hirose T. *et al*, *Eur J Biochem.* 1997 Oct **15**;249(2):481-8.

Description: CDK17/Cyclin Y complex of human **CDK17**, also known as PCK2, PCTAIRE2, GenBank Accession No. NM_002595, a.a. 1-523(end) with N-terminal GST-tag, MW=88 kDa, and human **Cyclin Y**, also known as CCNY, CBCP1, CCNX, and CFP1, GenBank Accession No. NM_145012, a.a. 2-341(end) with N-terminal GST-tag. MW=64 kDa. Co-expressed in a Baculovirus infected Sf9 cell expression system.

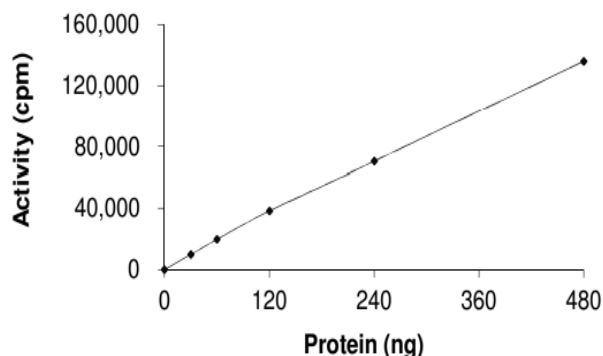
Specific Activity: ≥ 20 pmol/ $\mu\text{g}/\text{min}$

Assay Conditions: 5 mM MOPS, 2.5 mM β -glycerophosphate, 5 mM MgCl_2 , 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 50 ng/ml BSA, 2 mM ATP, 200 μg of MBP (or Histone H1) as substrate. Add [^{33}P]-ATP and incubate at 30°C for 15 minutes, then spot reaction on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



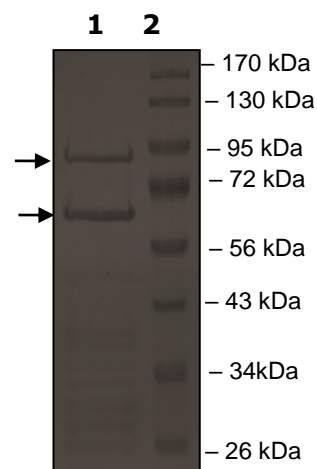
4-20% SDS-PAGE Coomassie staining

Lane 1:
CDK17/Cyclin Y

Lane 2:
Protein Marker

Purity: $\geq 70\%$

MW:
CDK17: 88 kDa
Cyclin Y: 64 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com