

Data Sheet

CDK9/CyclinT1, GST-tag

Human, recombinant, N-terminal GST-tag

Catalog #: 40307

Lot #: 140514 **Conc.:** 0.14 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 20% glycerol, and 1.6 mM glutathione

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Yang, Z., *et al.*, *Nat.* 2001; **414**: 317-322.
2. Baumli, S., *et al.*, *EMBO J.* 2008; **27(13)**: 1907-1918.

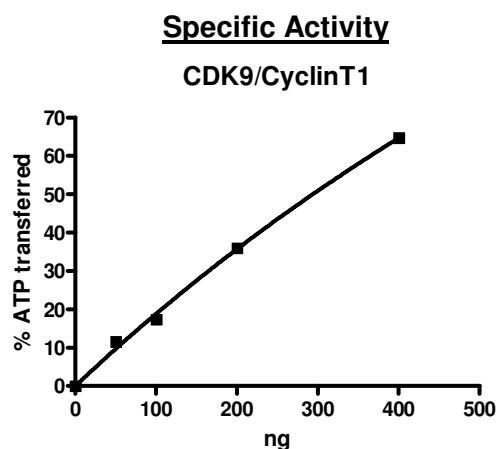
Description: Human cyclin-dependent kinase 9 (CDK9), also known as C-2K and cell division protein kinase 9, complexed with human cyclin T1 (CCNT1, CytT1), GenBank Accession Nos. NM_001261 (CDK9) and NM_001240 (CyclinT1), a.a. 2-372(end) (CDK9) and a.a. 2-726(end) (CyclinT1) each with N-terminal GST-tags, co-expressed in a baculovirus-infected Sf9 cell expression system. MW = 69 kDa (CDK9) and 107 kDa (CyclinT1).

Specific Activity: 26 pmol/min/ μg

Assay Conditions: Assay was done in a Kinase buffer using PDKtide (0.1 mg/ml) as a substrate with 10 μM ATP at 30°C for 45 min. The amount of ATP transferred from autophosphorylation as well as substrate phosphorylation was calculated using Kinase-Glo reagent (Promega).

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

Quality Assurance



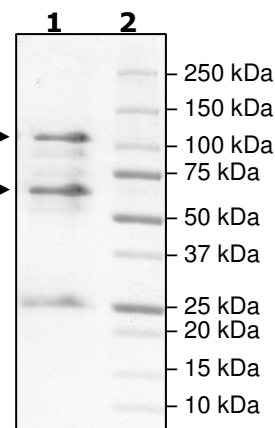
4-20% SDS-PAGE
Coomassie staining

Lane 1:
1.8 μg CDK9/CyclinT1

Lane 2:
Protein Marker

Purity: $\geq 66\%$

MW:
CyclinT1: 107 kDa
CDK9: 69 kDa



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com