

Data Sheet

CDK4/CyclinD3, GST-His-Tag

Human, recombinant, CDK4 N-terminal GST-His-Tag, CyclinD3 N-terminal GST-Tag
Catalog #: 40104
Lot #: 180710-3 **Conc.:** 0.11 mg/ml

Formulated in: 40mM Tris-HCl pH 8.0, 110mM NaCl, 2.2mM KCl, 20% Glycerol and 3mM DTT and 2 mM Glutathione

Stability: At least 6 months at -80°C

References:

1. Bullrich, F. *et al.*, *Cancer Res.* (1995); **55**:1199-1205.
2. Ye, X. *et al.*, *Proc. Nat. Acad. Sci.* (2001); **98**:1682-1686.

Description:

Complex of recombinant human CDK4, a.a. 1-303 (end), with N-Terminal GST-His-Tag, GenBank Accession No. NM_000075, and Cyclin D3, a.a 2-292(end), with N-terminal GST-tag, GenBank Accession No. NM_001760. Co-expressed in a Baculovirus-infected Sf9 cell expression system. CDK4 MW=64 kDa, Cyclin D3 MW=59 kDa.

Specific Activity: 10 pmole/min/ug.

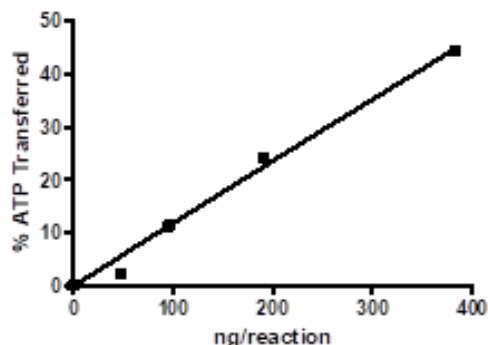
Assay Conditions: Assay was done in a Kinase buffer using Rb (0.02 mg/ml) as a substrate with 10 µM ATP at 30°C for 60 min. The amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).

Application:

Useful for three-dimensional structure studies.

Quality Assurance

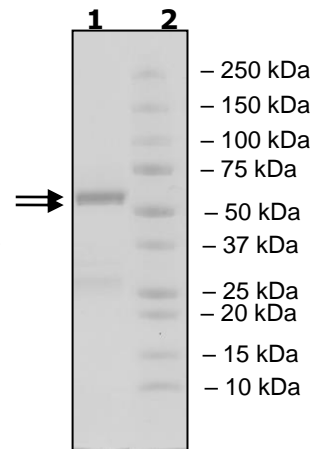
Specific Activity



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 µg Cdk4/CyclinD3
Lane 2:
Protein Marker

MW: 64/59kDa
Purity: 73%



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