

## Data Sheet

### CDK4(EE, T172A)/ Cyclin D1

Human, recombinant

**Catalog #:** 40094

**Lot #:** 100624

**Conc.:** 0.62 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
124 mM NaCl, 2.4 mM KCl, 225 mM  
imidazole, 10% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

**References:**

1. Amendola D. *et al.*, *Cell Prolif.* (2009) Feb; **42(1)**:94-109.
2. Yang L. *et al.*, *Biochem Cell Biol.* (2009) Dec; **87(6)**:933-42.

**Description:**

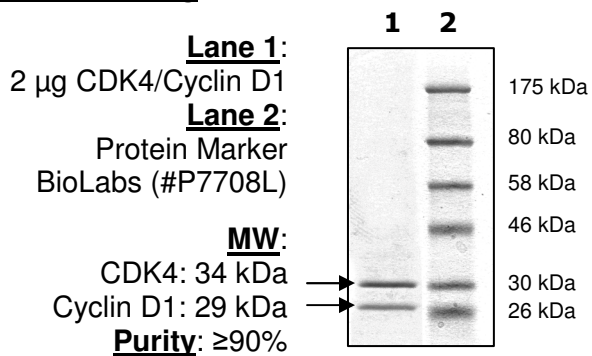
Complex of C-terminal His-tagged recombinant human CDK4 (GenBank Accession No. NM\_000075) (43G to 47G were replaced with EE and T172A), and Cyclin D1 (GenBank Accession No. NM\_053056) (a.a. 16 – 271). Co-expressed in a Baculovirus-infected Sf9 cell expression system. CDK4 MW=34 kDa, Cyclin D1 MW=29 kDa.

**Application:**

Useful for three-dimensional structure studies.

## Quality Assurance

4-20% SDS-PAGE  
Coomassie staining



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)