

## Data Sheet

### CDK3/CyclinE1

Human, recombinant, N-terminal GST tag

**Catalog #:** 40103

**Lot #:** 110520

**Conc.:** 0.1 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.1 mM EGTA, 0.1 mM EDTA, 0.1 mM PMSF, 25 % glycerol and 3 mM DTT.

**Stability:** >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.

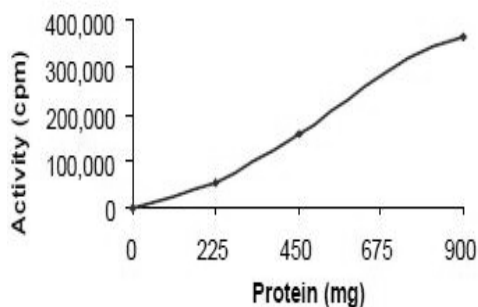
**Description:**

Complex of N-terminal GST-tagged recombinant full length human CDK3 (GenBank Accession No. NM\_001258), and Cyclin E1 (GenBank Accession No. NM\_001238). Co-expressed in a Baculovirus-infected Sf9 cell expression system. CDK3 MW=60 kDa, Cyclin E1 MW=73 kDa.

**Application:**

Useful for three-dimensional structure studies.

### Quality Assurance



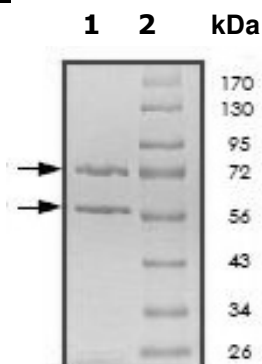
**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 CDK3/Cyclin E1

**Lane 2:**  
 Protein Marker

**MW:**  
 CDK3: 60 kDa  
 Cyclin E1: 73 kDa

**Purity:** ≥90%



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)