

## Data Sheet

### CDK2/CyclinE1, GST-Tag

Human, Recombinant, N-terminal GST-tag

**Catalog #:** 40102

**Lot #:** 171121      **Conc.:** 0.1 mg/ml

**Formulated in:** 50 mM Tris-HCl, 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

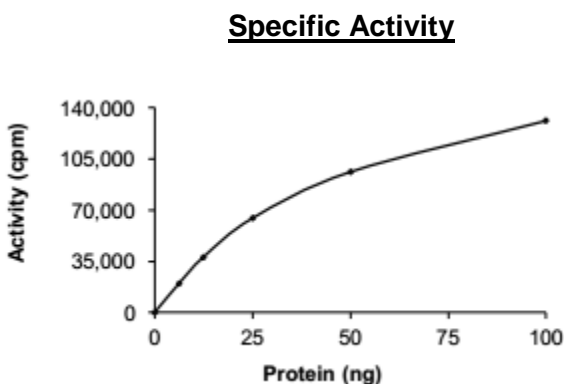
1. Caldon, C.E., *et al Cell Div.* 2010; **5(2)**.
2. Huang, H., *et al. Science* 2006; **314:294-297**.

**Description:** Human Dual CDK2/CyclinE1 protein complex also known as p33 and CCNE, respectively, GenBank Accession No. NM\_001798 and NM\_001238, both full length, with N-terminal GST-tags, MW=58 kDa (CDK2) and MW=73 kDa (CyclinE1), expressed in a baculovirus-infected Sf9 cell expression system.

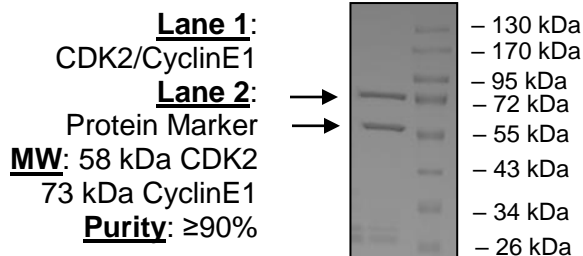
**Specific Activity:** 343 pmole/min/ $\mu\text{g}$   
**Assay Conditions:** Varying concentrations of CDK2/CyclinE1 were mixed with 5  $\mu\text{g}$  of Histone H1 protein substrate and 5  $\mu\text{l}$  of 250  $\mu\text{M}$  [ $^{33}\text{P}$ ]-ATP was added for a total volume of 25  $\mu\text{l}$ . Mixture was incubated for 15 min at  $30^{\circ}\text{C}$ . Reaction was stopped by placing 20  $\mu\text{l}$  of reaction onto phosphocellulose P81 paper. Paper was washed 3 times with 1% phosphoric acid solution for 10 min each and air dried. Radioactivity was measured from the P81 paper in the presence of scintillation fluid using a scintillation counter.

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

### 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)