

## Data Sheet

### CDK2(1-298)/CyclinA2

Human, recombinant, N-terminal His-GST tag  
 Catalog #: 41101  
 Lot #: 130325                      Conc.: 0.46 mg/ml

**Formulated in:** 20 mM NaH<sub>2</sub>PO<sub>4</sub>/Na<sub>2</sub>HPO<sub>4</sub>, pH 8.0, 400 mM NaCl, 20 mM Glutathione, 10% glycerol, and 2 mM DTT.

**Stability:** >6 months at -70°C. Avoid multiple freeze/thaw cycles.

**References:**

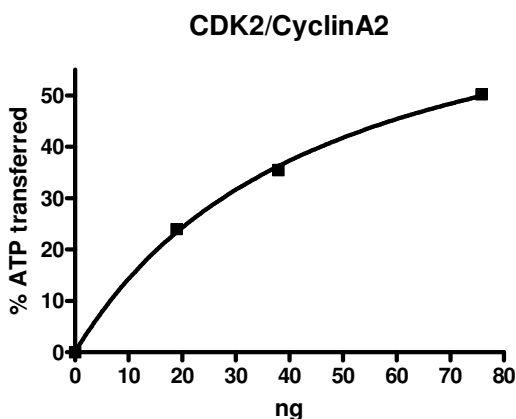
1. Brown NR., *et al. Cell Cycle*. 2007 Jun 1;**6(11)**:1350-9.
2. Hara KT., *et al. Dev Biol*. 2005 Oct 1;**286(1)**:102-13.

**Description:** Recombinant full-length human CDK2 a.a. 1-298 (end), MW = 34 kDa with no tag, and CyclinA2 a.a. 174-432 (end) with N-terminal His-GST tags, MW = 57 kDa. Both were co-expressed in a baculovirus Sf9 insect cell expression system. The CDK2 gene accession number is NM\_001798 and the CyclinA2 gene accession number is NM\_001237.

**Specific Activity:** 203 pmol/min/μg.  
**Assay Conditions:** Assay was done in a Kinase buffer using Histone H1 (Cell Signaling; 0.1 mg/ml at final concentration) as a substrate with 20 μM ATP at 30°C for 45 min. The amount of ATP transferred 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:**  
 3.48 μg CDK2/Cyclin A2

**Lane 2:**  
 Protein Marker

**MW:**  
 Cyclin A2= 57 kDa  
 CDK2= 34 kDa

**Purity:** ≥98%

