

# Certificate of Analysis



## CDK1

(cyclin dependent kinase 1)

Recombinant Human Protein Kinase

Synonyms: CDC2

Product No.: 0134-0000-1

Lot: 001

**Description:** Human CDK1, amino acids M<sub>1</sub>-M<sub>297</sub> (as in GenBank entry NM\_001786)\*, N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in Sf9 insect cells

\*Sequence may contain documented polymorphisms  
Detailed aa-sequence on request

**Product identity:** CDK1 was confirmed as CDK1 by specific Western Blotting

**Theoretical MW**<sub>GST-CDK1</sub>: 63,882 Da

**Expression:** Baculovirus infected Sf9 cells

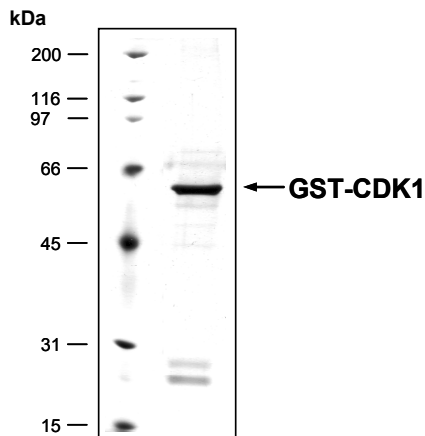
**Purification:** GST-Affinity Chromatography

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 7.5 mM reduced glutathione, 20% glycerol

**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

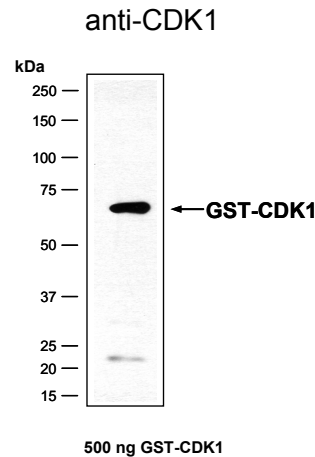
**Protein concentration:** 0.080 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

**Coomassie stain:**

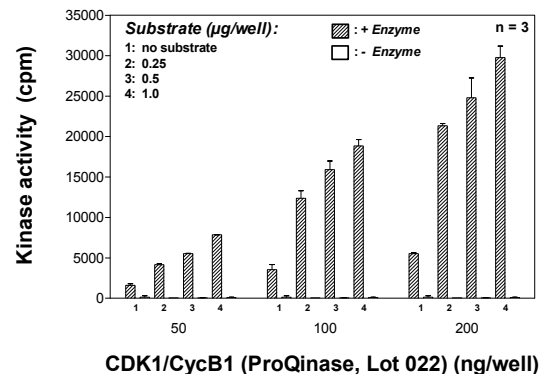
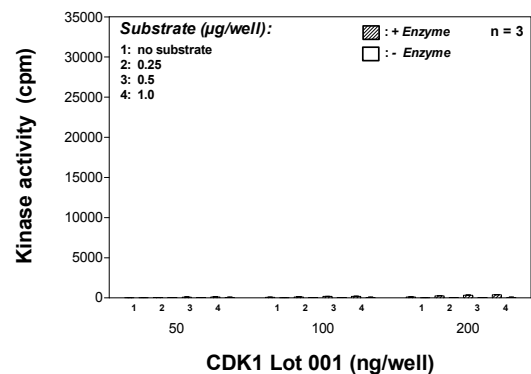


2.0 µg GST-CDK1

**Western blot analysis:**



**Kinase activity CDK1 vs active CDK1/CycB1:**



**Final assay concentrations:**

- 60 mM HEPES-NaOH, pH 7.5
- 3 mM MgCl<sub>2</sub>
- 3 mM MnCl<sub>2</sub>
- 3 µM Na-orthovanadate
- 1.2 mM DTT
- 50 µg / ml PEG<sub>20,000</sub>
- 1 µM ATP (610,000 cpm <sup>33</sup>P-ATP)
- Substrate (variable): Histone H1
- Recombinant CDK1 or CDK1/CycB1 (variable)

**Assay:** <sup>33</sup>PanQinase® Assay

This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.