

## Cdc25A

Cell division cycle 25A protein phosphatase  
human, recombinant, *E. coli*

Cat. No.	Amount
PR-348	20 $\mu$ g

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 25 mM HEPES pH 7.2, 40 mM NaCl and 2 mM DTT.

#### Assay buffer

50 mM Tris-HCl, pH 8.0, 50 mM NaCl, 0.1 mM EDTA.

#### Activity

1 nmol of Cdc25A will hydrolyze 10 nmol of pNPP per minute at 37°C at pH 8.0.

#### Molecular Weight

68 kDa

#### Purity

≥ 90% by SDS-PAGE

### Description

The dual specific Cdc25A (Cell division cycle) phosphatase regulates cell cycle progression, has oncogenic and anti-apoptotic activity, and is overexpressed in many human tumors.

### Selected References:

Fauman *et al.* (1998) Crystal structure of the catalytic domain of the human cell cycle control phosphatase, Cdc25A. *Cell* **93**:617.  
Gottlin *et al.* (1996) Kinetic analysis of the catalytic domain of human cdc25B. *J. Biol. Chem.* **271**:27445.