# **DATA SHEET**





### Cdc25A

Cell division cycle phosphatase human, recombinant, *E. coli* 

Cat. No.	Amount
PR-348	20 μg

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months Molecular Weight: 68 kDa Accession number: M81933 Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 20 mM Tris-HcL pH 8.0, 100 mM NaCl, 1 mM

EDTA and 1 mM DTT)

# 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. The protein does not have tags.

Assay buffer: 50 mM Tris-HCl, pH 8.0, 50 mM NaCl, 0.1 mM

## **Activity:**

1 nmol of Cdc25A will hydrolyze 10 nmol of pNPP per minute at  $37^{\circ}$ C at pH 8.0.

#### **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.