

CDC7/DBF4

(CDC7: Cell division cycle 7-related protein kinase, CDC7L1,
DBF4: Protein DBF4 homolog A, Activator of S phase kinase, Chiffon homolog A,
DBF4-type zinc finger-containing protein 1, ASK, DBF4A, ZDBF1)

- **CAT#:** CDC7/DBF4
- **Enzyme:** Human CDC7/DBF4
- **Substrate:** Peptide substrate, [KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYIADWC], 20 μ M
ATP 10 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	80
Ro 31-8220	37,000
GW 5074	1,200
H-89	>50,000
Rottlerin	>50,000

- **Scientific Information:** CDC7: O00311 (<http://www.expasy.org/uniprot/O00311>)
DBF4: Q9UBU7 (<http://www.expasy.org/uniprot/Q9UBU7>)

Reference compound IC50 for CDC7/DBF4

