

CDK9/Cyclin T1

(Cell division protein kinase 9, Cyclin-dependent kinase 9, Serine/threonine-protein kinase PITALRE, C-2K, Cell division cycle 2-like protein kinase 4, CDC2L4)

- **CAT#:** CDK9/Cyclin T1
- **Enzyme:** Human CDK9/Cyclin T1
- **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	14
Ro 31-8220	10
GW 5074	8,500
H-89	12,000

- **Scientific Information:** P50750 (<http://www.expasy.org/uniprot/P50750>)

Reference compound IC50 for CDK9/cyclinT1

