

---

## CDK9/CycT1

Product code 04-110

Full-length human CDK9 [1-372(end) amino acids of accession number NP\_001252.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-CyclinT1 [1-726(end) amino acids of accession number NP\_001231.2] using baculovirus expression system. GST-CDK9 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CDK9 substrate

ATP ( $\mu$ M) Km app / Bin : 9.4 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 5.2

IC50 at 1 mM ATP (nM) : 130

---