

---

---

## 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	: MSA
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

---

---