

---

## CDK7/CycH/MAT1

Product code 04-108

Full-length human CDK7 [1-346(end) amino acids of accession number NP\_001790.1] was co-expressed as N-terminal GST-fusion protein (66 kDa) with CyclinH [1-323(end) amino acids of accession number NP\_001230.1] and MAT1 [1-309(end) amino acids of accession number NP\_002422.1] using baculovirus expression system. GST-CDK7 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CTD3 peptide

ATP ( $\mu$ M) Km app / Bin : 32 / 50

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 17

IC50 at 1 mM ATP (nM) : 120

---