

---

---

# 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.

Assay platform : MSA  
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

---

---