

---

# CDK4/CycD1

Product code: 04-120

Full-length human CDK4 [1-303(end) amino acids of accession number NP\_000066.1] was co-expressed as N-terminal GST-fusion protein (61 kDa) with DYKDDDDK tagged CyclinD1(CCND1) [1-295(end) amino acids of accession number NP\_444284.1] using baculovirus expression system. GST-CDK4/CycD1 was purified by using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : DYRKtide-F

ATP ( $\mu$ M) Km app / Bin : 800 / 1000  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 72  
IC50 at 1 mM ATP (nM) : 72

---