

---

---

## CDK6/CycD3

Product code: 04-107

Full-length human CDK6 [1-326(end) amino acids of accession number NP\_001250.1] was co-expressed as N-terminal GST-fusion protein (64 kDa) with human GST-CyclinD3 [1-292(end) amino acids and S259A of accession number NP\_001751.1] using baculovirus expression system. GST-CDK6/CycD3 was purified by using glutathione sepharose chromatography and activated with His-CDK7/CycH/MAT1. Activated GST-CDK6/CycD3 was purified using glutathione sepharose chromatography.

Assay platform : MSA  
Substrate : DYRKtide-F  
ATP ( $\mu$ M) Km app / Bin : 330 / 300  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 58  
IC50 at 1 mM ATP (nM) : 110

---

---