

---

---

## CDK3/CycE1

Product code: 04-104

Full-length human CDK3 [1-305(end) amino acids of accession number NP\_001249.1] was co-expressed as N-terminal GST-fusion protein (62kDa) with CyclinE1 [1-410(end) amino acids of accession number NP\_001229.1] using baculovirus expression system.

Assay platform : MSA  
Substrate : Modified Histone H1  
ATP ( $\mu$ M) Km app / Bin : 1000 / 1000  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 3.4  
IC50 at 1 mM ATP (nM) : 3.4

---

---