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