
CDK4/CycD3

Product code 04-105

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 human GST-CyclinD3 [1-292(end) amino acids of accession number AAA51927.1] using baculovirus expression system. GST-CDK4/CycD3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : DYRKtide-F

ATP (μ M) Km app / Bin : 200 / 200

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 13

IC50 at 1 mM ATP (nM) : 52
