
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 and S259A of accession number NP_001751.1] using baculovirus expression system. GST-CDK4/CycD3 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
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

