
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 (μ M) Km app / Bin	: 330 / 300
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 58
IC50 at 1 mM ATP (nM)	: 110
