
CDK2/CycE1

Product code: 04-165

Full-length human CDK2 [1-298(end) amino acids of accession number NP_001789.2] was co-expressed as N-terminal GST-tagged protein (61 kDa) with CyclinE1 [1-410(end) amino acids of accession number NP_001229.1] using baculovirus expression system. GST-CDK2 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: Modified Histone H1
ATP (μ M) Km app / Bin	: 130 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.8
IC50 at 1 mM ATP (nM)	: 10
