
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.

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
