
CDK1(CDC2)/CycB1

Product code: 04-102

Full-length human CDK1 [1-297(end) amino acids of accession number NP_001777.1] was co-expressed as N-terminal GST-fusion protein (61 kDa) with CyclinB1 [1-433(end) amino acids of accession number NP_114172.1] using baculovirus expression system.

Assay platform : MSA
Substrate : Modified Histone H1
ATP (μ M) Km app / Bin : 34 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 3.3
IC50 at 1 mM ATP (nM) : 32
